Chromogranin A Recombinant Monoclonal Rabbit Antibody (CHGA/1731R)



Product Description

Chromogranin A is present in neuroendocrine cells throughout the body, including the neuroendocrine cells of the large and small intestine, adrenal medulla and pancreatic islets. It is an excellent marker for carcinoid tumors, pheochromocytomas, paragangliomas, and other neuroendocrine tumors. Co-expression of chromogranin A and neuron specific enolase (NSE) is common in neuroendocrine neoplasms. Reportedly, co-expression of certain keratins and chromogranin indicates neuroendocrine lineage. The presence of strong anti-chromogranin staining and absence of anti-keratin staining should raise the possibility of paraganglioma. The co-expression of chromogranin and NSE is typical of neuroendocrine neoplasms. Most pituitary adenomas and prolactinomas readily express chromogranin.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the CF® Dye Brochure for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1731, Anti-Chromogranin A (CHGA/1731R)

Product attributes

Call us: 800-304-5357

Antibody number #1731 Antibody reactivity (target) CHGA, Chromogranin A Antibody type Primary Host species Rabbit Clonality Recombinant Monoclonal Clone CHGA/1731R Isotype IgG, kappa Molecular weight 68-75 kDa Synonyms Beta-Granin; CGA; CHGA; Chromogranin A Parathyroid Secretory Protein 1; Fituitary Secretory Protein 1; SP-1; Vasostatin 1 or II Human gene symbol CHGA Entrez gene ID 1113 SwissProt P10645 Unigene 150793 Immunogen Recombinant human full-length Chromogranin A protein Verified antibody applications Verified antibody application Antibody target cellular localization Verified antibody application notes Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody application should be determined by user Positive control <th>Product attributes</th> <th></th>	Product attributes				
Antibody type Primary Host species Rabbit Clonality Recombinant Monoclonal Clone CHGA/1731R Isotype IgG, kappa Molecular weight 68-75 kDa Synonyms Beta-Granin; CGA; CHGA; Chromogranin A Parathyroid Secretory Protein 1; ER-37; Pancreastatin; Parathyroid Secretory Protein 1; ER-37; Pancreastatin; Parastatin; Parathyroid Secretory Protein 1; Pitultary Secretory Protein 1; Pitultary Secretory Protein 1; Phuman gene symbol CHGA Entrez gene ID 1113 SwissProt P10645 Unigene 150793 Immunogen Recombinant human full-length Chromogranin A protein Verified antibody applications Antibody target cellular localization Protein 1; Pitultary Secretory Protein 1; Pitultary Secret	Antibody number	#1731			
Host species Rabbit	Antibody reactivity (target)	CHGA, Chromogranin A			
Clonality Recombinant Monoclonal Clone CHGA/1731R Isotype IgG, kappa Molecular weight 68-75 kDa Synonyms Bela-Granin; CGA; CHGA; Chromogranin A Parathyroid Secretory Protein 1; En-3.7; Pancreastatin; Parastatin; Parathyroid Secretory Protein 1; Pituitary Secretory Protein 1; SP-I; Vasostatin 1 or II Human gene symbol CHGA Entrez gene ID 1113 SwissProt P10645 Unigene 150793 Immunogen Recombinant human full-length Chromogranin A protein Verified antibody applications HC (FFPE) (verified) Antibody target cellular localization Vesicular Species reactivity Human Antibody application notes Higher concentration may be required for direct detection with secondary antibody conjugates than for indirect detection with secondary antibody conjugates than for indirect detection with secondary antibody (ormalin): 0.25-0.5 u/gmll., Staining of formalin-liked tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 u/gmillion cells(0.1 mt., Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition	Antibody type	Primary			
Clone CHGA/1731R Isotype IgG, kappa Molecular weight 68-75 kDa Synonyms Beta-Granin; CGA; CHGA; Chromogranin A Parathyroid Secretory Protein 1; EP-37; Pancreastatin; Parathyroid Secretory Protein 1; EP-37; Pancreastatin; Parathyroid Secretory Protein 1; SP-1; Vasostatin 1 or II Human gene symbol CHGA Entrez gene ID 1113 SwissProt P10645 Unigene 150793 Immunogen Recombinant human full-length Chromogranin A protein Verified antibody applications HC (FFPE) (verified) Vesicular Vesicular Antibody application notes Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody (formalin): 0.25-0.5 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/mllion cells/0.1 mL, Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition Room temperature Storage Conditions Storage Conditions For research use only (RUO) Antibody/conjugate For Product may contain either bovine serum albumin (BSA) from PRS villow azide. Purified, BSA-free: 1 mg/mL in PBS villow azide. Purified, BSA-free: 1 mg/mL in PBS villow azide. Purified, BSA-free: 1 mg/mL in PBS villows a packed in Chinese hamster ovary cells. Inquire for the specific lot. Cell/tissue expression Neuroendocrine cells	Host species	Rabbit			
Isotype	Clonality	Recombinant Monoclonal			
Molecular weight Synonyms Beta-Granin; CGA; CHGA; Chromogranin A Parathyroid Secretory Protein 1; ER-37; Pancreastatin; Parastatin; Parathyroid Secretory Protein 1; Pituitary Secretory Protein 1; SP-I; Vasostatin I or II Human gene symbol CHGA Entrez gene ID 1113 SwissProt P10645 Unigene 150793 Immunogen Recombinant human full-length Chromogranin A protein Verified antibody applications Antibody target cellular localization Species reactivity Human Antibody application notes Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody conjugates than for indirect detection with secondary antibody (manunofluorescence: 1-2 ug/mL, Immunohistology (formalin): 0.25-0.5 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by conign at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition Room temperature Storage Conditions Storage Conditions Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate formulation PG0.05% BSA/0.05% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS without azide Validated in protein array Monospecific Antibody research areas Endocrinology, Neuroscience Product origin Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.	Clone	CHGA/1731R			
Beta-Granin; CGA; CHGA; Chromogranin A Parathyroid Secretory Protein 1; ER-37; Pancreastatin; Parastatin; Parathyroid Secretory Protein 1; Pituitary Secretory Protein 1; Prituitary Secretory Protein 1; Prituitary Secretory Protein 1; SP-1; Vasostatin 1 or II	Isotype	IgG, kappa			
Secretory Protein 1; ER-37; Pancreastatin; Parastatin; Parastatin; Parathyroid Secretory Protein 1; Pituitary Secretory Protein 1; SP-1; Vasostatin 1 or II Human gene symbol CHGA Entrez gene ID 1113 SwissProt P10645 Unigene 150793 Immunogen Recombinant human full-length Chromogranin A protein Verified antibody applications IHC (FFPE) (verified) Vesicular Vesicular Species reactivity Human Antibody application notes Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Immunohistology (formalin): 0.25-0.5 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition Room temperature Storage Conditions Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate formulation Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA,	Molecular weight	68-75 kDa			
Entrez gene ID SwissProt P10645 Unigene 150793 Immunogen Recombinant human full-length Chromogranin A protein Verified antibody applications Antibody target cellular localization Species reactivity Human Antibody application notes Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/ml. Immunohistology (formalin): 0.25-0.5 ug/ml., Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0 for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition Room temperature Storage Conditions Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate formulation For research use only (RUO) Antibody/conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.	Synonyms	Secretory Protein 1; ER-37; Pancreastatin; Parastatin; Parathyroid Secretory Protein 1; Pituitary Secretory Protein I;			
SwissProt P10645 Unigene 150793 Immunogen Recombinant human full-length Chromogranin A protein Verified antibody applications IHC (FFPE) (verified) Antibody target cellular localization Vesicular Species reactivity Human Antibody application notes Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescene: 1-2 ug/mL, Immunohistology (formalin): 0.25-0.5 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate formulation Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS	Human gene symbol	CHGA			
Unigene 150793 Immunogen Recombinant human full-length Chromogranin A protein Verified antibody applications IHC (FFPE) (verified) Antibody target cellular localization Vesicular Species reactivity Human Antibody application notes Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/ml., Immunohistology (formalin): 0.25-0.5 ug/ml., Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0 for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate formulation Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% azide, Purified; 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified; 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified; 0.2 mg/mL	Entrez gene ID	1113			
Immunogen Recombinant human full-length Chromogranin A protein Verified antibody applications IHC (FFPE) (verified) Antibody target cellular localization Vesicular Species reactivity Human Antibody application notes Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescene: 1-2 ug/ml, Immunohistology (formalin): 0.25-0.5 ug/ml, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate formulation Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.	SwissProt	P10645			
Verified antibody applications Antibody target cellular localization Species reactivity Human Antibody application notes Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/ml., Immunohistology (formalin): 0.25-0.5 ug/ml., Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition Room temperature Storage Conditions Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate formulation Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.	Unigene	150793			
applications Antibody target cellular localization Species reactivity Human Antibody application notes Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/mL, Immunohistology (formalin): 0.25-0.5 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition Room temperature Storage Conditions Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate formulation Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, PRSS/0.05% BSA/0.05% azide, PRSS/0.05% BSA/0.05% azide, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purifie	Immunogen	Recombinant human full-length Chromogranin A protein			
Species reactivity		IHC (FFPE) (verified)			
Antibody application notes Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/ml. Immunohistology (formalin): 0.25-0.5 ug/ml., Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition Room temperature Storage Conditions Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate Conjugates: 0.1 mg/mL in PBS/0.0% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% azide, Pu		Vesicular			
primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Immunoflistology (formalin): 0.25-0.5 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user Positive control PC12 cells. Adrenal gland, bowel, parathyroid, pancreas, or pheochromocytoma. Shipping condition Room temperature Storage Conditions Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified; 0	Species reactivity	Human			
pheochromocytoma. Shipping condition Room temperature Storage Conditions Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate formulation Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% BSA/0.05% BSA/ree: 1 mg/mL in PBS/0.05% BSA/ree: 1 mg/mL in PBS/0.05% azide, Purified on the PBS/0.05% BSA/ree: 1 mg/mL in PBS/0.05% azide, Purified on the PBS/0.05% BSA/ree: 1 mg/mL in PBS/0.05% azide, Purified on the PBS/0.05% BSA/ree: 1 mg/mL in PBS/0.05% azide, Purified on the PBS/0.05% BSA/ree: 1 mg/mL in PBS/0.05% azide, Purified on the PBS/0.05% BSA/ree: 1 mg/mL in PBS/0.05% azide, Purified on the PBS/0.05% BSA/0.05% azide, Purified on PBS/0.05% BSA/0.05% BSA/0.05% azide, Purified on PBS/0.05% BSA/0.05% azide, Purified on PBS/0.05% BSA/0.05% BSA/0.05% azide, Purified on PBS/0.05% BSA/0.05% BSA/0.0	Antibody application notes	primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Immunohistology (formalin): 0.25-0.5 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal			
Storage Conditions	Positive control				
Note: store BSA-free antibodies at -10 to -35 °C Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate formulation Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide Validated in protein array Monospecific Antibody research areas Endocrinology, Neuroscience Product origin Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. Cell/tissue expression Neuroendocrine cells	Shipping condition	Room temperature			
stored as recommended Regulatory status For research use only (RUO) Antibody/conjugate formulation Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide Validated in protein array Monospecific Antibody research areas Endocrinology, Neuroscience Product origin Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. Cell/tissue expression Neuroendocrine cells	Storage Conditions				
Antibody/conjugate formulation Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% BSA/0.05% BSA/0.05% BSA/free: 1 mg/mL in PBS without azide Validated in protein array Monospecific Antibody research areas Endocrinology, Neuroscience Product origin Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. Cell/tissue expression Neuroendocrine cells	Shelf life				
tormulation conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% BSA/0.05% BSA/0.05% BSA/0.05% BSA/free: 1 mg/mL in PBS/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS/0.05% azide, Purified, DSA-free: 1 mg/mL in PBS/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS/0.05% azide, Purified, P	Regulatory status	For research use only (RUO)			
Antibody research areas Endocrinology, Neuroscience Product origin Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. Cell/tissue expression Neuroendocrine cells		conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL ir			
Product origin Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. Cell/tissue expression Neuroendocrine cells	Validated in protein array	Monospecific			
bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. Cell/tissue expression Neuroendocrine cells	Antibody research areas	Endocrinology, Neuroscience			
·	Product origin	bovine serum (Bos taurus), or recombinant BSA produced in			
Tumor expression Neuroendocrine cancer, Pituitary cancer	·				
	Tumor expression	Neuroendocrine cancer, Pituitary cancer			

Email: btinfo@biotium.com

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on November 3, 2025 at 04:10:23 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/recombinant-rabbit-monoclonal-anti-chromogranin-a-chga-chga1731r/