## HER-2 / CD340 Monoclonal Mouse Antibody (ERBB2/2452)



## **Product Description**

This antibody recognizes a protein of 185 kDa, which is identified as c-erbB-2/HER-2/neu. Its epitope is localized in the extracellular domain. C-erbB-2/HER-2 is a member of the EGFR family. This MAb is specific and shows minimal cross-reaction with other members of the EGFR-family. Receptors of this family are located on the plasma membrane and consist of an extracellular ligand-binding domain that is connected to a large intracellular domain by a single transmembrane sequence. c-erbB-2/HER-2 protein is over-expressed in a variety of carcinomas especially those of breast and ovary.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 2452, Anti-HER-2 / c-ErbB-2 / Neu / CD340 (ERBB2/2452)

Call us: 800-304-5357

## Product attributes

Product attributes				
Antibody number	#2452			
Antibody reactivity (target)	c-ErbB-2, CD340, HER-2, Neu			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	ERBB2/2452			
Isotype	IgG2a, kappa			
Molecular weight	185 kDa			
Synonyms	p185, CD340, Verb b2 Erythroblastic Leukemia Viral Oncogene Homolog 2, Neuro/Glioblastoma Derived Oncogene Homolog			
Human gene symbol	ERBB2			
Entrez gene ID	2064			
SwissProt	P04626			
Unigene	446352			
Immunogen	Recombinant human HER-2 protein fragment (around aa 311-462) (exact sequence is proprietary)			
Verified antibody applications	Flow (verified), IF (verified), IHC (FFPE) (verified)			
Antibody target cellular localization	Plasma membrane			
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, Immunohistology (formalin): 1-2 ug/mL for 30 minutes at RT, Staining of formalin-fixed tissues requires			
	boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user			
Positive control	minutes followed by cooling at RT for 20 minutes, Optimal			
Positive control Shipping condition	minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user			
	minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user SKBR-3 Cells or Breast Cancers.			
Shipping condition	minules followed by cooling at RT for 20 minules, Optimal dilution for a specific application should be determined by user SKBR-3 Cells or Breast Cancers.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Shipping condition Storage Conditions	minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user SKBR-3 Cells or Breast Cancers.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when			
Shipping condition Storage Conditions Shelf life	minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user SKBR-3 Cells or Breast Cancers.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user SKBR-3 Cells or Breast Cancers.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  For research use only (RUO)  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; 0.2 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 in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user SKBR-3 Cells or Breast Cancers.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  For research use only (RUO)  Conjugates: 0.1 mg/mL in PBS/0.05% 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			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation Validated in protein array	minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user SKBR-3 Cells or Breast Cancers.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  For research use only (RUO)  Conjugates: 0.1 mg/mL in PBS/0.05% 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  Monospecific			

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,	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 30, 2025 at 02:12:16 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/her-2-c-erbb-2-neu-cd340-monoclonal-mouse-antibody-erbb2-2452/">https://biotium.com/product/her-2-c-erbb-2-neu-cd340-monoclonal-mouse-antibody-erbb2-2452/</a>