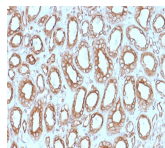


Calnexin Monoclonal Mouse Antibody (CANX/1543)



Product Description

Calnexin is a type-I integral membrane protein of the endoplasmic reticulum (ER). It is a calcium-binding chaperone protein that interacts transiently with newly synthesized N-linked glycoproteins, facilitating protein folding and assembly. It may also play a central role in the quality control of protein folding by retaining incorrectly folded protein subunits within the ER for degradation.

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.

Product attributes

Antibody number	#1543
Antibody reactivity (target)	Calnexin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CANX/1543
Isotype	IgG1, kappa
Molecular weight	67 kDa (predicted); 80-90 kDa (observed)
Synonyms	Calnexin; CANX; IP90; Major histocompatibility complex class I antigen-binding protein p88
Human gene symbol	CANX
Entrez gene ID	821
SwissProt	P27824
Unigene	567968
Immunogen	Recombinant N-terminal fragment of human Calnexin protein (around aa 1-300) (exact sequence is proprietary)
Antibody target cellular localization	Endoplasmic reticulum
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Species reactivity	Human
Positive control	HeLa, MCF-7 or U2OS cells. Kidney or small intestine.
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, Purified, BSA-free: 1 mg/mL in PBS without azide
Validated in protein array	Monospecific
Antibody research areas	Chaperones, Immunology