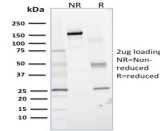


# Lactoylglutathione Lyase Monoclonal Mouse Antibody (CPTC-GLO1-1)



## Product Description

GLO1 is an enzyme involved in the detoxification of methylglyoxal, a byproduct of glycolysis. GLO1 expression has been demonstrated by several studies to be upregulated in various human malignant tumors, including metastatic melanoma and lung carcinoma, and thus is a target for pharmaceutical development.

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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

## Product attributes

Antibody number	#2258
Antibody reactivity (target)	Lactoylglutathione Lyase
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CPTC-GLO1-1
Isotype	IgG1, kappa
Molecular weight	24-26 kDa
Synonyms	Aldoketomutase; epididymis secretory protein Li 74; Glo1; Glx I; glyoxalase 1 complex; glyoxalase 1 regulatory; glyoxalase 1 structural; glyoxalase domain containing 1; Glyoxalase I; glyoxylase 1; Ketone-aldehyde mutase; lactoyl glutathione lyase; Lactoylglutathione lyase; Methylglyoxalase; S-D-lactoylglutathione methylglyoxal lyase
Human gene symbol	GLO1
Entrez gene ID	2739
SwissProt	Q04760
Unigene	268849
Immunogen	Recombinant human full-length GLO1 protein
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Cytoplasmic
Species reactivity	Human
Positive control	A431, HepG2, MOLT and Raji cell lysates, Ovarian carcinoma or prostate carcinoma.
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
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Tumor expression	Lung cancer, Melanoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
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, or LI-COR Bioscience.