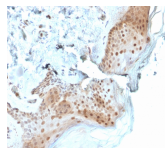


8-Oxoguanine DNA Glycosylase Monoclonal Mouse Antibody (CPTC-OGG1-1)



Product Description

8-Oxoguanine DNA Glycosylase (OGG1), otherwise known as N-Glycosylase/DNA lyase, is the enzyme responsible for the excision of 8-oxoguanine (8-oxoG), a mutagenic base by-product which occurs as a result of exposure to reactive oxygen. The action of this enzyme includes lyase activity for chain cleavage. Accumulation of 8-oxoG increases the occurrence of A:T to C:G or G:C to T:A transversion mutations, because 8-oxoG forms a stable basepair with adenine as well as with cytosine. The OGG1 gene encodes eight isoforms (OGG1A-C, OGG2A-E) which result from alternative splicing of a single messenger RNA. The OGG1A splice variant is the most prevalent form and localizes to the nucleus, whereas the OGG2A splice variant is targeted to the mitochondria. Defects in OGG1 may be a cause of renal cell carcinoma. 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 2251, 8-Oxoguanine DNA Glycosylase Monoclonal Mouse Antibody (CPTC-OGG1-1)

Product attributes

Antibody number	#2251
Antibody reactivity (target)	N-Glycosylase/DNA lyase
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CPTC-OGG1-1
Isotype	IgG2c, kappa
Molecular weight	39 kDa
Synonyms	8-oxoguanine DNA glycosylase; HMMH; HOGG1; MUTM; OGH1
Human gene symbol	OGG1
Entrez gene ID	4968
SwissProt	O15527
Unigene	380271
Immunogen	Recombinant full-length human OGG1 protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Mitochondria, Nucleus
Species reactivity	Human
Positive control	HeLa or Jurkat cells. Kidney, skin or lymph node.
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended

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
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, LI-COR Bioscience.