

Mouse Anti-PEG mAb

[Anti-Polyethylene glycol monoclonal antibody]

Catalog Number:	57000
Concentration:	1mg/ml
Origin of Antibody:	Mouse
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Mouse IgM in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	PEG conjugated to KLH
Purification:	Affinity chromatography.
Specificity:	The antibody has high affinity for the PEG backbone. The antibody binds to the variety of PEG such as PEG40K, PEG20K, PEG5K, PEG12, PEGylated drugs and PEG conjugates.
Applications:	ELISA Capturer: 5-10 ug/mL, ELISA Detector: 0.1-1.0 ug/mL, Western Blot: 0.1-1.0 ug/mL, Immunohistochemistry: 10-15 ug/mL
Note:	Some reagents such as Tween-20, Triton X-100 and NP-40 interfere with the PEG antibody interaction due to structural similarity to PEG. Antibody solution should not contain such reagents in the assay of PEGylated molecules.