

TP53BP1 Monoclonal Antibody [1B9]

[Catalog No. A-6118-100]

Description

Mouse monoclonal antibody raised against a partial recombinant TP53BP1, clone 1B9.

Size

100 µg

Purification

Protein A purified

Immunogen

TP53BP1 (NP_005648, 1766 a.a. ~ 1875 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG1 Kappa

Formulation

In 1x PBS, pH 7.2

Storage

Aliquot and store at -20°C or lower. Avoid repeated freezing and thawing. Multiple freeze/thaw cycles may result in decreased performance. Non-hazardous. No MSDS necessary.

Handling Recommendations

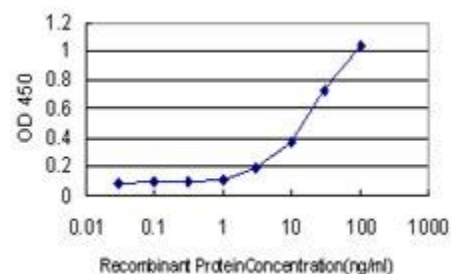
For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Applications & Suggested Dilutions*

ELISA: 0.25-1 µg/ml

**The end user is responsible for determining optimal working dilutions.*

Data



Detection limit for recombinant GST tagged TP53BP1 is approximately 0.1 ng/ml as a capture antibody.

This product is for research purposes only. Not intended for use in diagnostic procedures.