

GSTP1 Monoclonal Antibody [2G6-F6]

[Catalog No. A-6193-100]

Description

Mouse monoclonal antibody raised against a full length recombinant GSTP1, clone 2G6-F6.

Size

100 µg

Purification

Protein A purified

Immunogen

GSTP1 (AAH10915, 1 a.a. ~ 211 a.a) full length 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*

Western Blot: 0.5-1 µg/ml

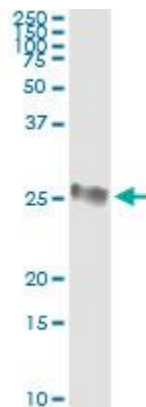
Immunoprecipitation: 5 µg/ml

ELISA: 0.25-1 µg/ml

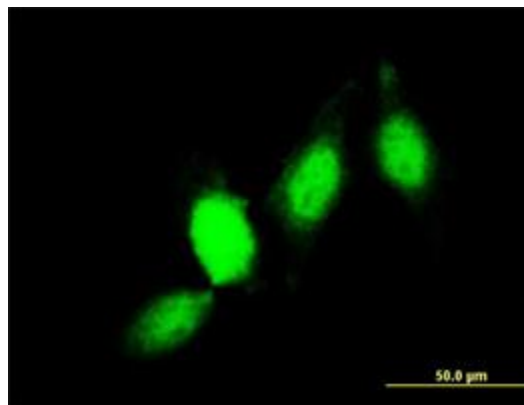
Immunofluorescence: 5 µg/ml

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

Data



Immunoprecipitation of GSTP1 transfected lysate using an anti-GSTP1 monoclonal antibody and Protein A Magnetic Bead, and immunoblotted with GSTP1 MaxPab rabbit polyclonal antibody.



Immunofluorescence of monoclonal antibody to GSTP1 on HeLa cell. [Antibody Concentration: 10 µg/ml]

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