

Gstp1 Polyclonal Antibody

(Catalog A56198)

Background

Conjugation of reduced glutathione to a wide number of exogenous and endogenous hydrophobic electrophiles. Can metabolize 1-chloro-2,4-dinitrobenzene. Regulates negatively CDK5 activity via p25/p35 translocation to prevent neurodegeneration.

Description

Gstp1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P19157

Purification

>95%, Protein G purified

Immunogen

Recombinant Mouse Glutathione S-transferase P 1 protein (2-210AA)

Storage

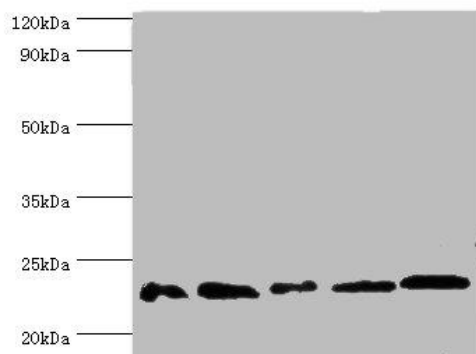
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

GST YF-YF, GST class-pi, GST-piB, Preadipocyte growth factor, Gstpib, Gstp1

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: Gstp1 Polyclonal Antibody at 2ug/ml

Lane 1: mouse kidney tissue

Lane 2: jurkat whole cell lysate

Lane 3: k562 whole cell lysate

Lane 4: 293T whole cell lysate

Lane 5: Hela whole cell lysate

Secondary

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 23kDa

Observed band size: 23kDa