
GFAP Polyclonal Antibody, HRP Conjugated

(Catalog # A52934)

Background

GFAP, a class-III intermediate filament, is a cell-specific marker that, during the development of the central nervous system, distinguishes astrocytes from other glial cells.

Description

GFAP Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P14136

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Glial fibrillary acidic protein (292-432AA)

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

GFAP

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
