
DBI Polyclonal Antibody, Biotin Conjugated

(Catalog #A55163)

Background

Binds medium- and long-chain acyl-CoA esters with very high affinity and may function as an intracellular carrier of acyl-CoA esters. It is also able to displace diazepam from the benzodiazepine (BZD) recognition site located on the GABA type A receptor. It is therefore possible that this protein also acts as a neuropeptide to modulate the action of the GABA receptor.

Description

DBI Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P07108

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Acyl-CoA-binding protein (2-105AA)

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

ACBP, Diazepam-binding inhibitor, DBI, Endozepine, EP, DBI

Application

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