

Prod. no. Cyp-11

Product Description

RABBIT ANTI-FISH CYP1A PEPTIDE

Polyclonal antibody, CP-226

biosense
LABORATORIES

www.biosense.com

Description

This product consists of affinity-purified rabbit polyclonal antibodies against peptides 190-204 and 282-296 of rainbow trout cytochrome P450 1A (CYP1A).

The antibody is dissolved in 0.1 M Na-phosphate, 0.15 M NaCl, 1% BSA, 0.1% Na-azide, pH 7.4.

Immunogen source

Synthetic peptides corresponding to amino acid sequence 190-204 and 282-296 of rainbow trout (*Oncorhynchus mykiss*) CYP1A, the latter coupled to the carrier protein keyhole limpet hemocyanin (KLH).

Specificity

The monoclonal antibody Cyp-11 cross-reacts strongly to CYP1A in liver samples from a wide variety of species, including rainbow trout (*Oncorhynchus mykiss*), Atlantic salmon (*Salmo salar*), Atlantic cod (*Gadus morhua*), common carp (*Cyprinus carpio*), flounder (*Platichthys flesus*), sheepshead minnow (*Cyprinodon variegatus*) and gilthead seabream (*Sparus aurata*). The degree of cross-reactivity differs between species and with the methods employed.

Application

The polyclonal antibody Cyp-11 works well in both ELISA and western blot for a number of fish species. Since assay conditions vary, the optimum dilution should be determined for each particular application.

Normal dilution range:

ELISA: 1:100 - 1:2000

Western blot: 1:100 - 1:2000

Storage

In the dark at 4°C.

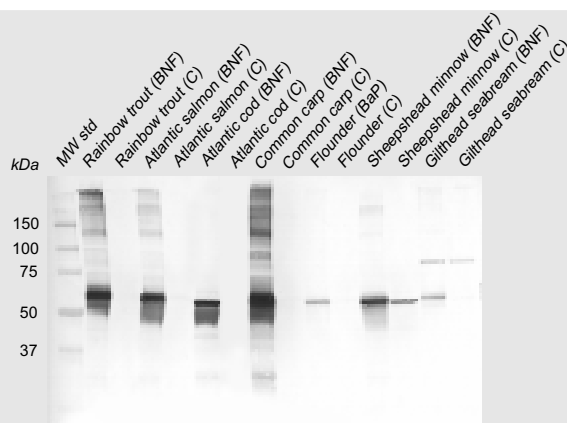
Western blot

Samples: Liver microsomes (10 µg protein/ml) from β-naphtoflavone (BNF) or benzo(a)pyrene (BaP) induced fish, or control fish (C).

Primary antibody: Cyp-11 diluted 1:500

Secondary antibody: Peroxidase labelled Goat anti-rabbit IgG, diluted 1:3000

Development: Peroxidase color development reagent (4-chloro-1-naphthol/H₂O₂)



ELISA

Coating: Liver microsomes (10 µg protein per well) from β-naphtoflavone (BNF) or benzo(a)pyrene (BaP) induced fish, or control fish (C).

Primary antibody: Cyp-11

Secondary antibody: Peroxidase labelled Goat anti-rabbit IgG, diluted 1:3000

Development: Peroxidase substrate (OPD/H₂O₂)

