

Mouse anti-brown bullhead Vtg Monoclonal antibody, ND-1D12

General description

This product consists of affinity-purified mouse monoclonal antibodies against vitellogenin from brown bullhead (*Ameiurus nebulosus*).

Immunogen source: Vitellogenin purified from brown bullhead (*Ameiurus nebulosus*), sheepshead minnow (*Cyprinodon variegatus*), bowfin (*Amia calva*) and spotted gar (*Lepisosteus oculatus*).

Subclass/isotype: IgG₁-kappa.

Specificity

The monoclonal antibody ND-1D12 binds with high affinity to vitellogenin from brown bullhead (*Ameiurus nebulosus*). The antibody does not cross-react with vitellogenin from sheepshead minnow (*Cyprinodon variegatus*), bowfin (*Amia calva*), spotted gar (*Lepisosteus oculatus*), killifish (*Fundulus heteroclitus*), tilapia (*Oreochromis niloticus*), halibut (*Hippoglossus hippoglossus*), rainbow trout (*Oncorhynchus mykiss*), carp (*Cyprino carpio*) or zebrafish (*Danio rerio*).

The degree of cross-reactivity may differ between species and with the methods employed.

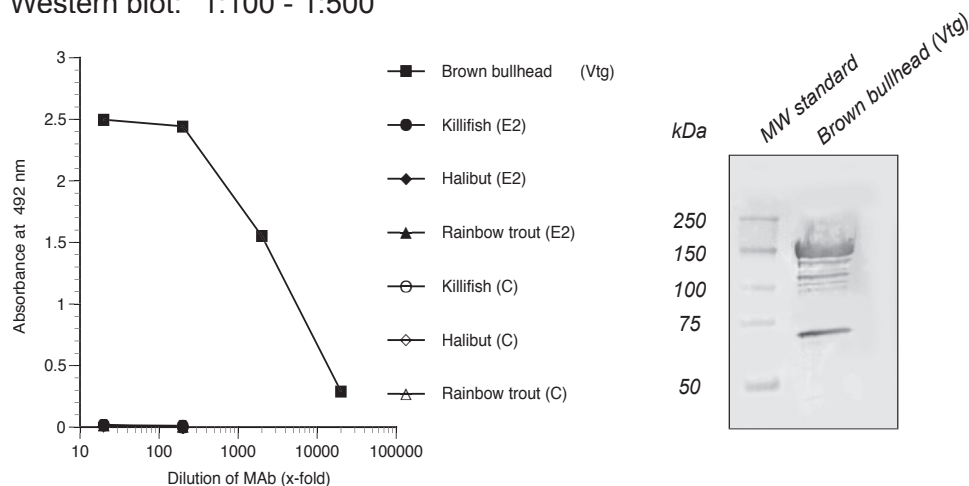
Applications

The monoclonal antibody ND-1D12 works well in both ELISA and western blot. Since assay conditions vary, the optimum dilution should be determined for each particular application. *Note*: The antibody is dissolved in a buffer containing BSA, and is therefore not recommended to use for coating.

Normal dilution range:

ELISA: 1:100 - 1:1000

Western blot: 1:100 - 1:500



ELISA

Coating: Plasma (pl; diluted 1:1000) from 17 β -estradiol treated (E2) or control (C) fish. Purified vitellogenin (Vtg; 2 μ g/ml).

Primary antibody: ND-1D12

Western blot

Samples: Purified vitellogenin (Vtg; 1 μ g/well).

Primary antibody: ND-1D12 diluted 1:200