

Mouse anti-carp Vtg Monoclonal antibody, ND-2D3

General description

This product consists of affinity-purified mouse monoclonal antibodies against vitellogenin from common carp (*Cyprinus carpio*).

Immunogen source: Vitellogenin purified from plasma of 17 β -estradiol treated striped bass (*Morone saxatilis*), brown bullhead (*Ameiurus nebulosus*) and common carp (*Cyprinus carpio*).

Subclass/isotype: IgG₁-kappa.

Specificity

The monoclonal antibody ND-2D3 binds with high affinity to vitellogenin from common carp (*Cyprinus carpio*). The antibody also cross-reacts with vitellogenin from fathead minnow (*Pimephales promelas*), zebrafish (*Danio rerio*), goldfish (*Carassius auratus*), mullet (*Mugil* sp.), pinfish (*Lagodon rhomboides*) and sucker (*Catostomidae* sp.). The antibody does not react with vitellogenin from striped bass (*Morone saxatilis*) or brown bullhead (*Ictalurus nebulosus*). The degree of cross-reactivity differs between species and with the methods employed.

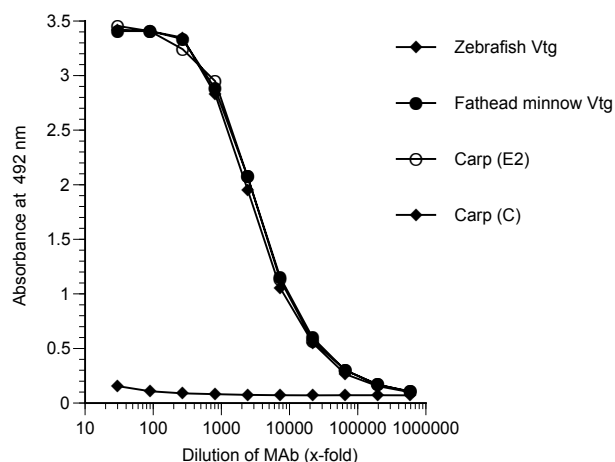
Applications

The monoclonal antibody ND-2D3 works well in both ELISA and western blot for a number of species. 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

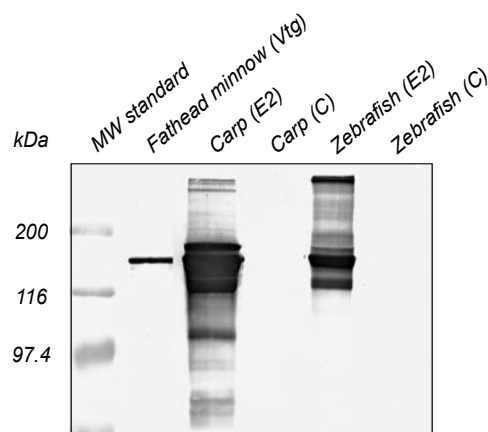
Western blot: 1:500 - 1:1000



ELISA

Coating: Plasma (diluted 1:1000) from 17 β -estradiol treated (E2) or control (C) fish. Purified vitellogenin (Vtg; 500 ng/well).

Primary antibody: ND-2D3



Western blot

Samples: Plasma (1:200, 10 μ l/well) from 17 β -estradiol treated (E2) or control (C) fish. Purified vitellogenin (Vtg; 5 μ g/well).

Primary antibody: ND-2D3 diluted 1:500