

## Mouse anti-striped bass Vtg Monoclonal antibody, ND-1C8

### General description

This product consists of affinity-purified mouse monoclonal antibodies against vitellogenin from striped bass (*Morone saxatilis*).

Immunogen source: Vitellogenin purified from plasma of 17 $\beta$ -estradiol treated striped bass (*Morone saxatilis*).

Subclass/isotype: IgG<sub>1</sub>-kappa.

### Specificity

The monoclonal antibody ND-1C8 binds with high affinity to vitellogenin from striped bass (*Morone saxatilis*) and rainbow trout (*Oncorhynchus mykiss*). The antibody also cross-reacts with vitellogenin from a variety of different species including largemouth bass (*Micropterus salmonides*), Atlantic cod (*Gadus morhua*), tilapia (*Oreochromis niloticus*), killifish (*Fundulus heteroclitus*) and sheepshead minnow (*Cyprinodon variegatus*). The degree of cross-reactivity differs between species and with the methods employed.

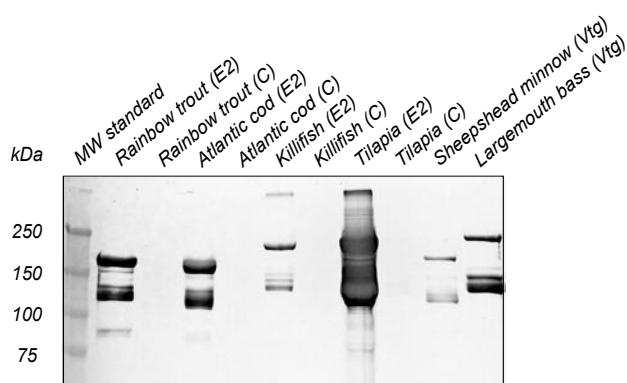
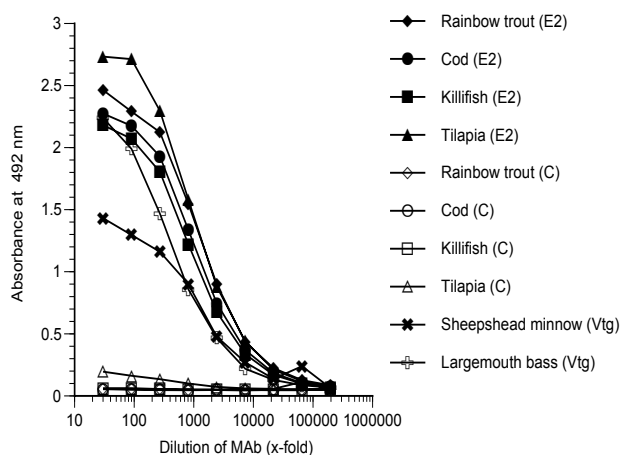
### Applications

The monoclonal antibody ND-1C8 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 $\beta$ -estradiol treated (E2) or control (C) fish. Purified vitellogenin (Vtg; 500 ng/well)

Primary antibody: ND-1C8

### Western blot

Samples: Plasma (1:200, 15  $\mu$ l/well) from 17 $\beta$ -estradiol treated (E2) or control (C) fish. Purified vitellogenin (Vtg; 1  $\mu$ g/well)

Primary antibody: ND-1C8 diluted 1:3000