

Mouse anti-killifish Vtg Monoclonal antibody, ND-5F8

General description

This product consists of affinity-purified mouse monoclonal antibodies against vitellogenin from killifish (*Fundulus heteroclitus*) and white perch (*Morone americanus*).

Immunogen source: Vitellogenin purified from plasma of 17 β -estradiol treated killifish (*Fundulus heteroclitus*), white perch (*Morone americanus*) and bowfin (*Amia calva*). Egg yolk from swordfish (*Xiphias gladius*).

Subclass/isotype: IgG₁-kappa.

Specificity

The monoclonal antibody ND-5F8 binds with high affinity to vitellogenin from killifish (*Fundulus heteroclitus*) and white perch (*Morone americanus*). The antibody also cross-reacts with vitellogenin from a variety of different species including common carp (*Cyprinus carpio*), wrasse (*Ctenolabrus rupestris*), tilapia (*Oreochromis niloticus*), halibut (*Hippoglossus hippoglossus*) and sheepshead minnow (*Cyprinodon variegatus*). The antibody does not cross-react with vitellogenin from bowfin (*Amia calva*) or with egg yolk from swordfish (*Xiphias gladius*). The degree of cross-reactivity differs between species and with the methods employed.

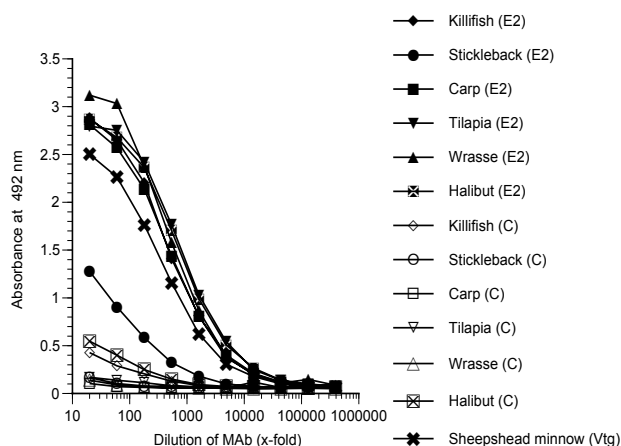
Applications

The monoclonal antibody ND-5F8 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

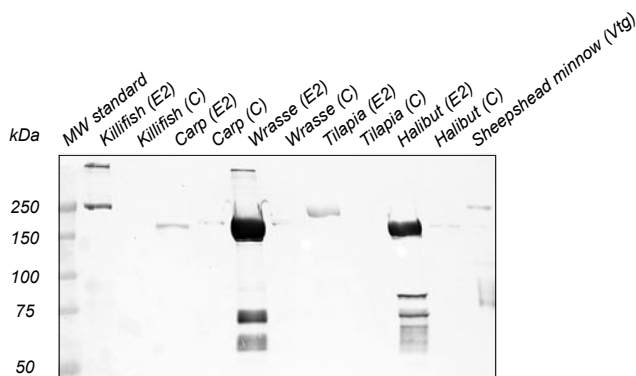
Western blot: 1:100 - 1:500



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-5F8



Western blot

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

Primary antibody: ND-5F8 diluted 1:100