

Mouse anti-sheepshead minnow Vtg Monoclonal antibody, ND-5C9

General description

This product consists of affinity-purified mouse monoclonal antibodies against vitellogenin from sheepshead minnow (*Cyprinodon variegatus*).

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

Subclass/isotype: IgG₁-kappa.

Specificity

The monoclonal antibody ND-5C9 binds with high affinity to vitellogenin in plasma samples from sheepshead minnow (*Cyprinodon variegatus*). The antibody also cross-reacts with vitellogenin from halibut (*Hippoglossus hippoglossus*) and flounder (*Platichthys flesus*). The antibody does not cross-react with vitellogenin from bowfin (*Amia calva*), brown bullhead (*Ameiurus nebulosus*) or spotted gar (*Lepisosteus oculatus*). The degree of cross-reactivity differs between species and with the methods employed.

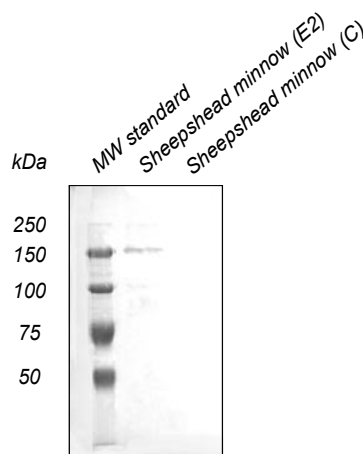
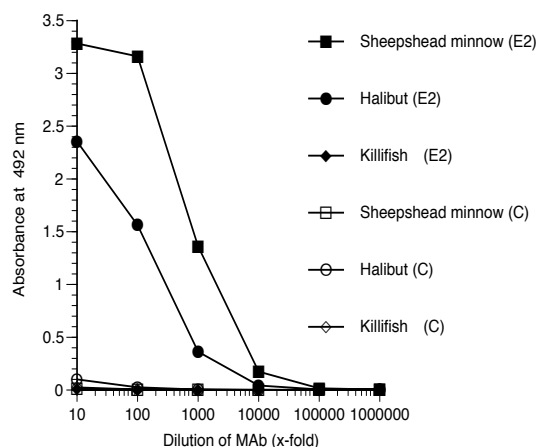
Applications

The monoclonal antibody ND-5C9 works well in ELISA, and can also be used in 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: 1:50 - 1:500



ELISA

Coating: Plasma (diluted 1:1000) from 17 β -estradiol treated (E2) or control (C) fish.

Primary antibody: ND-5C9

Western blot

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

Primary antibody: ND-5C9 diluted 1:50