

Mouse anti-salmon Vtg Monoclonal antibody, BN-5

General description

This product consists of affinity-purified mouse monoclonal antibodies against vitellogenin from Atlantic salmon (*Salmo salar*).

Immunogen source: Vitellogenin purified from plasma of 17 β -estradiol treated Atlantic salmon (*Salmo salar*).

Subclass/isotype: IgG₁-kappa.

Specificity

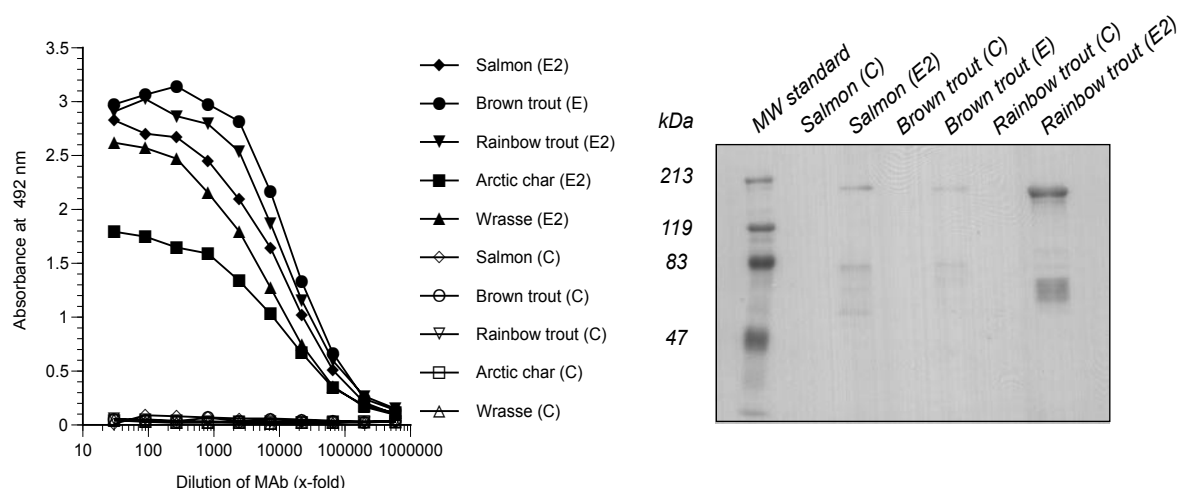
The monoclonal antibody BN-5 binds with high affinity to vitellogenin in plasma samples from Atlantic salmon (*Salmo salar*). The antibody cross-reacts with vitellogenin in plasma samples from brown trout (*Salmo trutta*), rainbow trout (*Oncorhynchus mykiss*), Arctic char (*Salvelinus alpinus*) and wrasse (*Ctenolabrus rupestris*). The degree of cross-reactivity differs between species and with the methods employed.

Applications

The monoclonal antibody BN-5 works well in ELISA for a number of species. The antibody may also work in western blot for some fish species, but the affinity for vitellogenin is much lower when this method is used. 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 - 1:2000



ELISA

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

Primary antibody: BN-5

Western blot

Samples: Plasma (1:200, 7 μ l/well) from 17 β -estradiol treated (E2), sexually mature female (E) or control (C) fish.

Primary antibody: BN-5 diluted 1:100