

Mouse anti-salmon Zrp Monoclonal antibody, MN-7F2

General description

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

Immunogen source: Purified eggshell proteins (*zona radiata* proteins and small amounts of *zona pellucida* proteins) from Atlantic salmon (*Salmo salar*).

Subclass/isotype: IgG₁-kappa.

Specificity

The monoclonal antibody MN-7F2 binds with high affinity to *zona radiata* proteins (Zrp) in plasma samples from Atlantic salmon (*Salmo salar*), reacting primarily with the γ monomer. The antibody also cross-reacts with Zrp from rainbow trout (*Oncorhynchus mykiss*). The degree of cross-reactivity differs between species and with the methods employed.

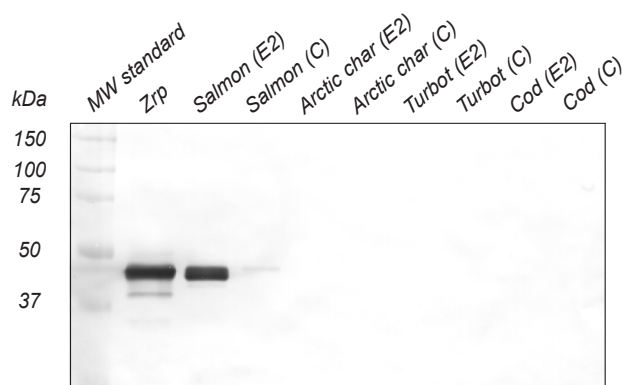
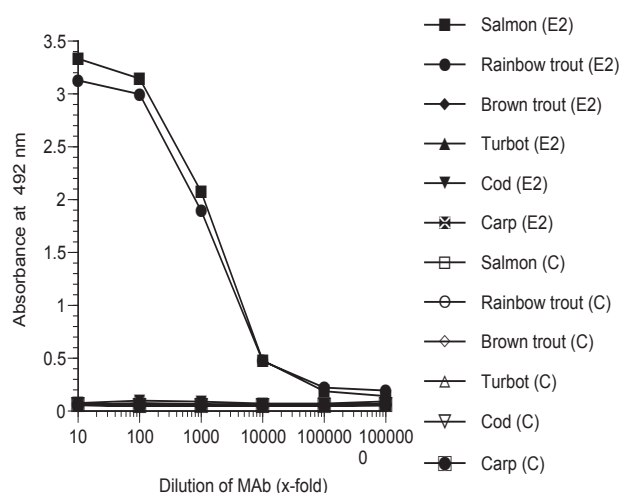
Applications

The monoclonal antibody MN-7F2 works well in both ELISA and 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 - 1:5000

Western blot: 1:100 - 1:1000



ELISA

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

Primary antibody: MN-7F2

Western blot

Samples: Purified α , β and γ Zrp, plasma (1:200, 10 μ l/well) from 17 β -estradiol treated (E2) or control (C) fish.

Primary antibody: MN-7F2 diluted 1:500