

## Mouse anti-salmon Zrp Monoclonal antibody, MN-8C4

### 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: IgG<sub>1</sub>-kappa.

### Specificity

The monoclonal antibody MN-8C4 binds with high affinity to *zona radiata* proteins (Zrp) in plasma samples from Atlantic salmon (*Salmo salar*), reacting primarily with the  $\alpha$  and  $\beta$  monomers. The antibody also cross-reacts with Zrp from rainbow trout (*Oncorhynchus mykiss*), Arctic charr (*Salvelinus alpinus*), turbot (*Scophthalmus maximus*) and cod (*Gadus morhua*). The degree of cross-reactivity differs between species and with the methods employed.

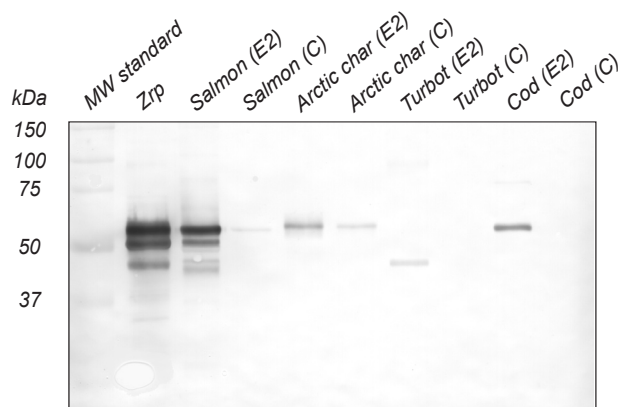
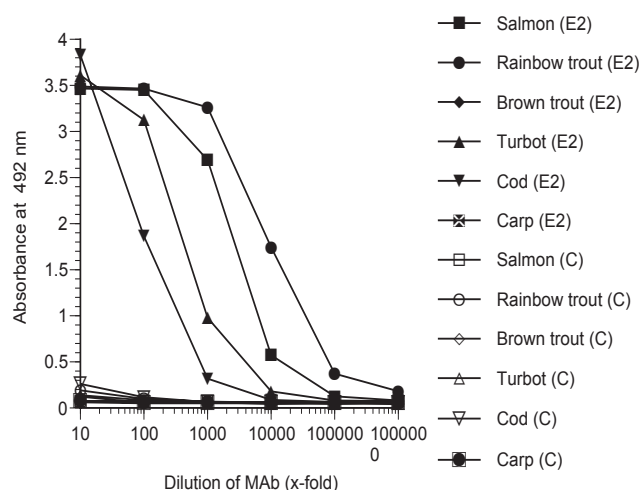
### Applications

The monoclonal antibody MN-8C4 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 - 1:5000

Western blot: 1:100 - 1:1000



### ELISA

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

Primary antibody: MN-8C4

### Western blot

Samples: Purified  $\alpha$ ,  $\beta$  and  $\gamma$  Zrp, plasma (1:200, 10  $\mu$ l/well) from 17 $\beta$ -estradiol treated (E2) or control (C) fish.

Primary antibody: MN-8C4 diluted 1:500