

Rabbit anti-sea bream Vtg Polyclonal antibody, PO-2

General description

This product consists of rabbit polyclonal antibodies (affinity-purified IgG fraction) against vitellogenin from sea bream (*Sparus aurata*).

Immunogen source: Vitellogenin purified from plasma of 17 β -estradiol treated sea bream (*Sparus aurata*).

Specificity

The polyclonal antibody PO-2 binds to vitellogenin in plasma samples from sea bream (*Sparus aurata*). The antibody also cross-reacts with vitellogenin from Atlantic salmon (*Salmo salar*), brown trout (*Salmo trutta*), rainbow trout (*Oncorhynchus mykiss*), turbot (*Scophthalmus maximus*), flounder (*Platichthys flesus*), Atlantic cod (*Gadus morhua*), wrasse (*Ctenolabrus rupestris*), tilapia (*Oreochromis niloticus*) and carp (*Cyprinus carpio*). The degree of cross-reactivity differs between species and with the methods employed.

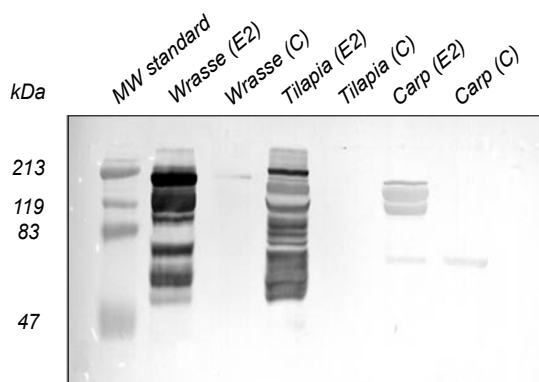
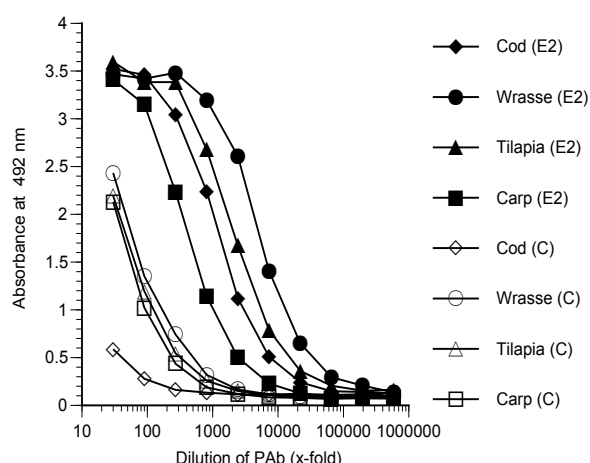
Applications

The polyclonal antibody PO-2 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: 1:200 - 1:3000

Western blot: 1:200 - 1:3000



ELISA

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

Primary antibody: PO-2

Western blot

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

Primary antibody: PO-2 diluted 1:3000