

Rabbit anti-cod Vtg Polyclonal antibody, CS-1

General description

This product consists of rabbit polyclonal antibodies (affinity-purified IgG fraction) against vitellogenin from Atlantic cod (*Gadus morhua*).

Immunogen source: Vitellogenin purified from plasma of 17 β -estradiol treated Atlantic cod (*Gadus morhua*).

Specificity

The polyclonal antibody CS-1 binds with high affinity to vitellogenin in plasma samples from Atlantic cod (*Gadus morhua*). The antibody also cross-reacts with vitellogenin from turbot (*Scophthalmus maximus*), rainbow trout (*Oncorhynchus mykiss*), common carp (*Cyprinus carpio*) and tilapia (*Oreochromis niloticus*). The degree of cross-reactivity differs between species and with the methods employed.

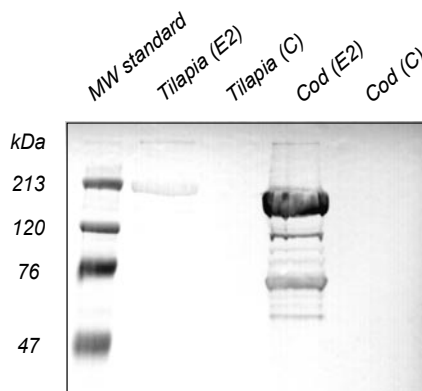
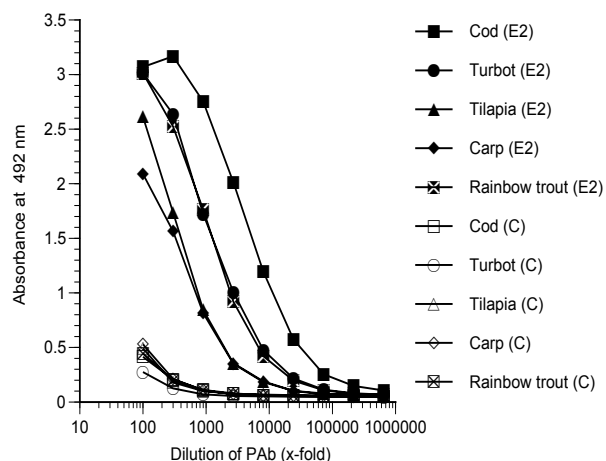
Applications

The polyclonal antibody CS-1 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 - 1:2000

Western blot: 1:200 - 1:3000



ELISA

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

Primary antibody: CS-1

Western blot

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

Primary antibody: CS-1 diluted 1:3000