

Rabbit anti-salmon Vtg Polyclonal antibody, AA-1

General description

This product consists of rabbit polyclonal antibodies (affinity-purified IgG fraction) against vitellogenin from Atlantic salmon (*Salmo salar*).

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

Specificity

The polyclonal antibody AA-1 binds with high affinity to vitellogenin in plasma samples from Atlantic salmon (*Salmo salar*). The antibody also cross-reacts with vitellogenin from other salmonids (*Salmoniformes*), flatfish (*Pleuronectiformes*), and Atlantic cod (*Gadus morhua*). The degree of cross-reactivity differs between species and with the methods employed.

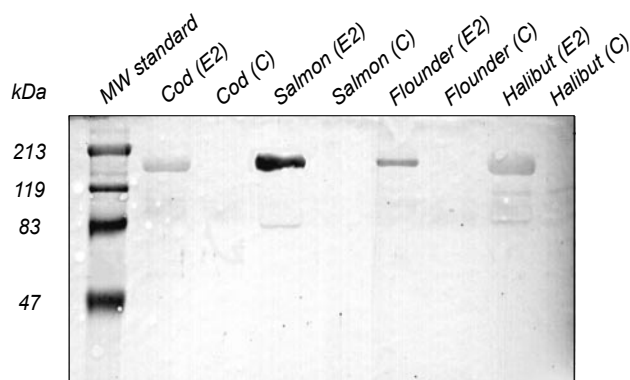
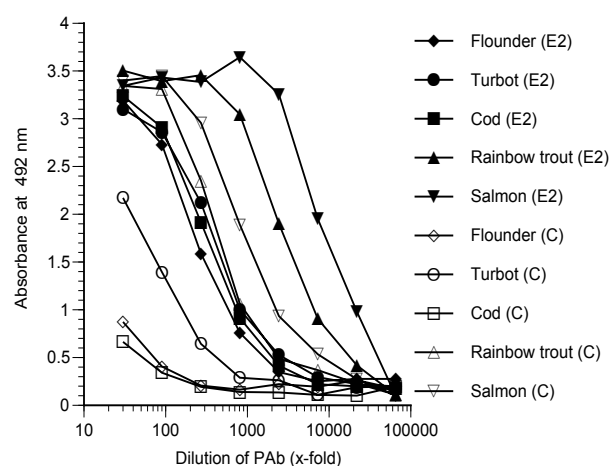
Applications

The polyclonal antibody AA-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: 1:50 - 1:3000

Western blot: 1:50 - 1:3000



ELISA

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

Primary antibody: AA-1

Western blot

Samples: Plasma (1:100, 10 μ l/well) from 17 β -estradiol treated (E2) or control (C) fish.

Primary antibody: AA-1 diluted 1:2000