

Mouse anti-zebrafish Vtg Monoclonal antibody, JE-10D4

General description

This product consists of affinity-purified mouse monoclonal antibodies against vitellogenin from zebrafish (*Danio rerio*).

Immunogen source: Vitellogenin purified from whole body homogenate of 17 β -estradiol treated zebrafish (*Danio rerio*).

Subclass/isotype: IgG₁-kappa.

Specificity

The monoclonal antibody JE-10D4 binds with high affinity to vitellogenin in whole body homogenate samples from zebrafish (*Danio rerio*). The antibody also cross-reacts with vitellogenin from other cyprinid species, including carp (*Cyprinus carpio*), fathead minnow (*Pimephales promelas*) and roach (*Rutilus rutilus*). The degree of cross-reactivity differs between species and with the methods employed.

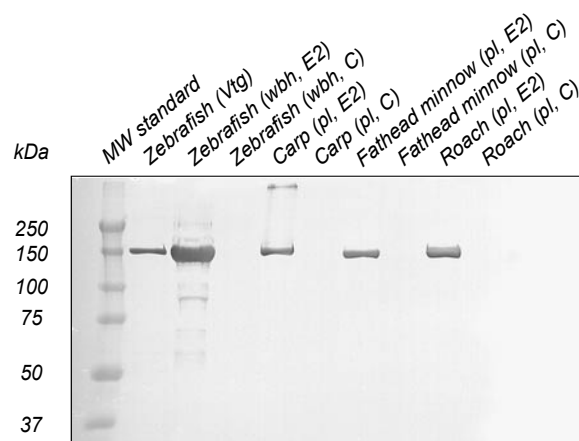
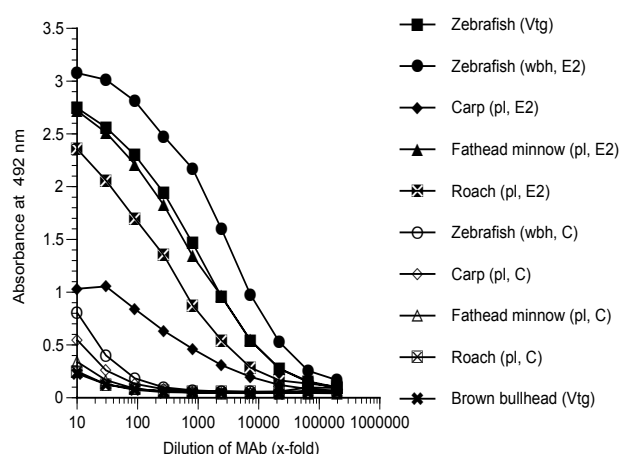
Applications

The monoclonal antibody JE-10D4 works well in both ELISA and western blot for a number of cyprinid species. The antibody also works well in immunohistochemistry. 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 - 1:1000

Western blot: 1:100 - 1:1000



ELISA

Coating: Whole body homogenate (wbh; 10 μ g protein/ml) or plasma (pl; diluted 1:1000) from 17 β -estradiol treated (E2) or control (C) fish. Purified vitellogenin (Vtg; 2 μ g/ml).

Primary antibody: JE-10D4

Western blot

Samples: Whole body homogenate (wbh; 10 μ g protein/well) or plasma (pl; 1:200, 10 μ l/well) from 17 β -estradiol treated (E2) or control (C) fish. Purified vitellogenin (Vtg; 1 μ g/well).

Primary antibody: JE-10D4 diluted 1:500