

## Mouse anti-Gulf sturgeon Vtg Monoclonal antibody, ND-1H2

### General description

This product consists of affinity-purified mouse monoclonal antibodies against vitellogenin from Gulf sturgeon (*Acipenser oxyrinchus desotoi*).

Immunogen source: Vitellogenin purified from plasma of 17 $\beta$ -estradiol treated Gulf sturgeon (*Acipenser oxyrinchus desotoi*).

Subclass/isotype: IgG<sub>1</sub>-kappa.

### Specificity

The monoclonal antibody ND-1H2 binds strongly to vitellogenin from Gulf sturgeon (*Acipenser oxyrinchus desotoi*). The antibody also cross-reacts with vitellogenin from lake sturgeon (*Acipenser fulvescens*). The degree of cross-reactivity may differ between species and with the methods employed.

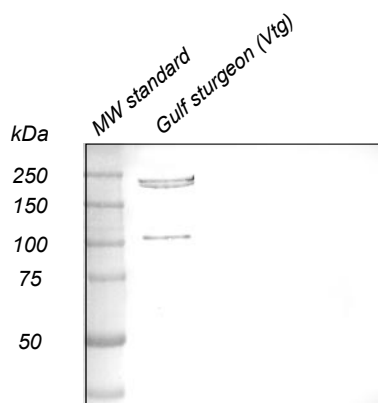
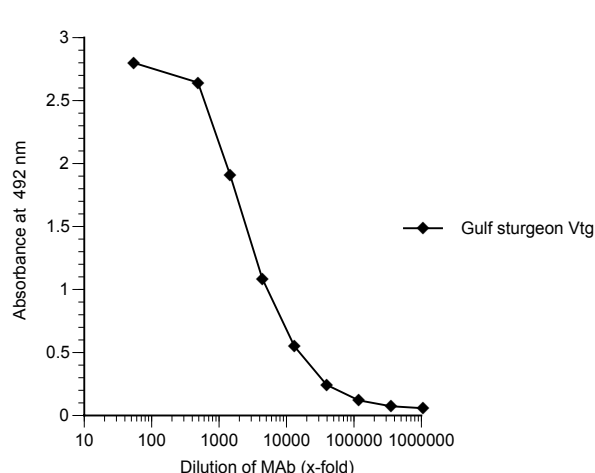
### Applications

The monoclonal antibody ND-1H2 binds to vitellogenin from Gulf sturgeon in both ELISA and western blot. 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 - 1:500



### ELISA

Coating: Purified vitellogenin (Vtg; 2  $\mu$ g/ml)

Primary antibody: ND-1H2

### Western blot

Sample: Purified vitellogenin (Vtg; 1  $\mu$ g/well)

Primary antibody: ND-1H2 diluted 1:100