

## Mouse anti-alligator Vtg Monoclonal antibody, ND-1E8

### General description

This product consists of affinity-purified mouse monoclonal antibodies against vitellogenin from alligator.

Immunogen source: Vitellogenin from American alligator (*Alligator mississippiensis*).

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

### Specificity

The monoclonal antibody ND-1E8 binds with high affinity to vitellogenin from American alligator (*Alligator mississippiensis*). The degree of cross-reactivity may differ between species and with the methods employed.

### Applications

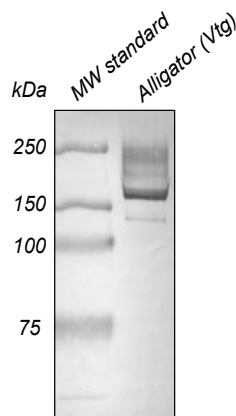
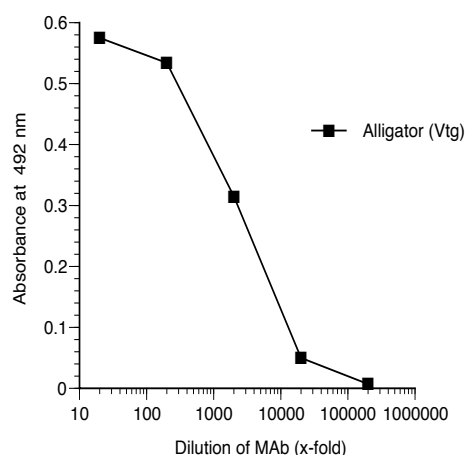
The monoclonal antibody ND-1E8 works well in western blot, and can also be used in ELISA. Since assay conditions vary, the optimum dilution should be determined for each particular application.

*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: Purified Vtg (2 µg/ml)

Primary antibody: ND-1E8

### Western blot

Sample: Purified Vtg (5 µg/well).

Primary antibody: ND-1E8 diluted 1:100