

Mouse anti-bird Vtg Monoclonal antibody, ND-3C3

General description

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

Immunogen source: Mixture of Vitellogenin from Japanese quail (*Coturnix japonica*), common tern (*Sterna hirundo*) and barn swallow (*Hirundo rustica*).

Subclass/isotype: IgG_{2b}-kappa.

Specificity

The monoclonal antibody ND-3C3 binds with high affinity to vitellogenin from Japanese quail (*Coturnix japonica*), common tern (*Sterna hirundo*) and barn swallow (*Hirundo rustica*). The antibody also cross-reacts with vitellogenin from other bird species, including American kestrel (*Falco sparverius*), zebra finch (*Taeniopygia guttata*) and domestic chicken (*Gallus gallus*). The degree of cross-reactivity differs between species and with the methods employed.

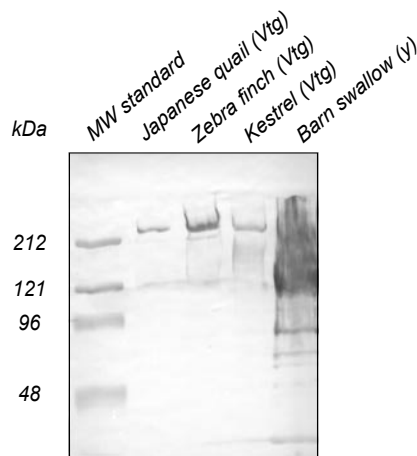
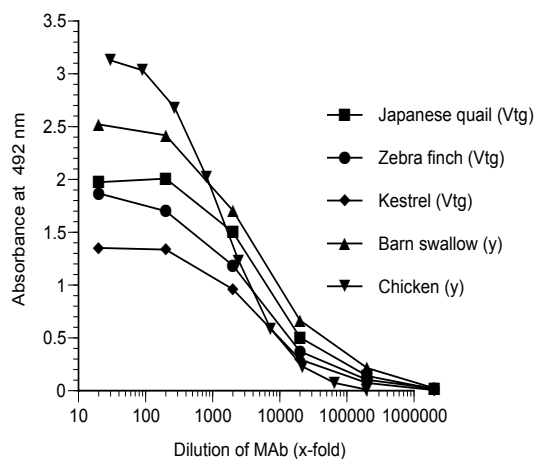
Applications

The monoclonal antibody ND-3C3 works well in both ELISA and western blot for a number of bird 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: 1:100 - 1:1000

Western blot: 1:50 - 1:500



ELISA

Coating: Purified Vtg (2 µg/ml) or egg yolk (y; diluted 1:500).

Primary antibody: ND-3C3

Western blot

Samples: Purified Vtg (5 µg/well) or egg yolk (y; diluted 1:500, 10 µl/well). Please note that the Vtg is somewhat degraded.

Primary antibody: ND-3C3 diluted 1:50