

Prod. no. Zrp-12

## Product Description

# RABBIT ANTI-SALMON ZONA RADIATA PROTEINS

Polyclonal IgG-fraction, O-146

**biosense**  
LABORATORIES

[www.biosense.no](http://www.biosense.no)

## Description

This product contains polyclonal rabbit antibodies (purified IgG fraction) against zona radiata proteins (Zrp) from Atlantic salmon. The antibody is dissolved in 0.1M Na-phosphate, 0.15M NaCl, 1% BSA, 0.1% Na-azide, pH 7.4.

## Immunogen source

Purified eggshell proteins (zona radiata proteins and small amounts of zona pellucida proteins) from Atlantic salmon (*Salmo salar*).

## Specificity

The polyclonal antibody O-146 binds with high affinity to the three zona radiata proteins  $\alpha$ ,  $\beta$  and  $\gamma$  in plasma samples from Atlantic salmon. The antibody cross-reacts well with Zrp in plasma samples from rainbow trout (*Oncorhynchus mykiss*) and Arctic char (*Salvelinus alpinus*), Atlantic cod (*Gadus morhua*), turbot (*Scophthalmus maximus*), flounder (*Platichthys flesus*) and halibut (*Hippoglossus hippoglossus*). The degree of cross-reactivity may differ between the different Zrp proteins and with the different methods used. Crossreactivity with Zona pellucida proteins may be observed in fish ovaries or eggs.

## Application

The polyclonal antibody Zrp-12 is well suited for ELISA and immunoblotting. Since assay conditions vary, the optimum dilution should be determined for each particular application.

Usual dilution range:

1:100 - 1: 1000 (*Salmoniformes*)

1:50 - 1: 200 (Other species)

## Storage

In the dark at 4 °C.

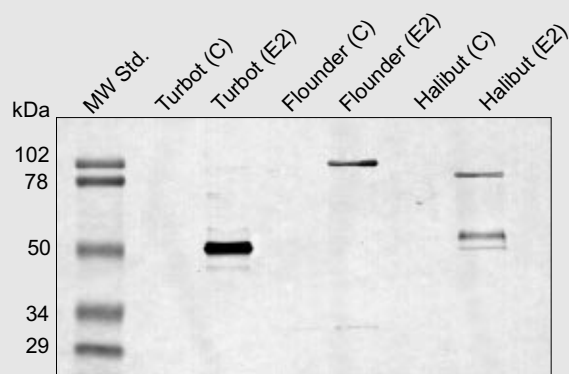
## Western blot

**Samples:** Plasma (1:100 dilution, 10  $\mu$ l) from juvenile fish induced with 17 $\beta$ -estradiol (E2) or control fish (C), non-induced juvenile or male fish.

**Primary antibody:** Zrp-12 diluted 1:100

**Secondary antibody:** Peroxidase labelled Goat anti-rabbit IgG, diluted 1:3000

**Development:** Peroxidase color development reagent (4-chloro-1-naphthol/H<sub>2</sub>O<sub>2</sub>)



## ELISA

**Coating:** Plasma (diluted 1:1000) from fish induced with 17 $\beta$ -estradiol (E2). Control fish (C) were non-induced juvenile or male fish.

**Primary antibody:** Zrp-12

**Secondary antibody:** Peroxidase labelled Goat anti-rabbit IgG, diluted 1:3000

**Development:** Peroxidase substrate (OPD/H<sub>2</sub>O<sub>2</sub>)

