

**Background:** Gonadotropin-releasing hormone (GnRH) is considered to stimulate LH secretion from the pituitary. This antibody is raised against an GnRH analog (GnRHa)

**Immunogen:** GnRHa conjugated to BSA.

**Application information**

Potent GnRH agonist. Induce stimulation of LH secretion from the pituitary and spawning activity in fish.

**ELISA:** 1:5000

In this immunoassay the anti-GnRHa and sample/standard are added simultaneously within GnRHa coated wells. After incubation, the anti-GnRHa bound to the coated GnRHa is detected with HRP-labelled secondary antibodies and color development achieved by addition of TMB reagent.

**Cross reactivity:**

N/A

**Product information:**

**Clonality:** Polyclonal

**Raised in:** Hen

**Purity:** Total IgY

**Antibody form:** Liquid

**Storage instructions:**

4°C - **DO NOT FREEZE!**

*Preparation contains sodium azide (0,02%) as inhibitor of bacterial growth.*

**Please,** remember to spin tubes briefly prior to opening them to avoid any losses that might occur from liquid or lyophilized material adhering to the cap or the sides of the tubes.

**Developed and produced by:**

AgriSera AB  
Box 57  
911 21 Vännäs  
Sweden

***For research use only, not for diagnostic or therapeutic use.***

[www.biosense.com](http://www.biosense.com)

Biosense Laboratories AS, Thormøhlensgt. 55, N-5008 Bergen, Norway  
Phone: +47 55543966, Fax: +47 55543771, e-mail: [biosense@biosense.no](mailto:biosense@biosense.no)