

# Product Information

logo

## Estrogen receptor $\alpha$ (F region) Monoclonal Antibody

Cat No: G01009 - 100  $\mu$ L

### General Data

<b>Shipping:</b>	wet ice
<b>Stability:</b>	Store at -20degrees; shelf life 3 year maximum after production Avoid repeat freezing and thawing cycles. Does not contain any preservative
<b>Formulation:</b>	Liquid Ascite, does not contain any preservative therefore avoid repeat freezing and thawing cycles
<b>Host:</b>	Mouse
<b>Antigen:</b>	synthetic peptide corresponding to amino acids 578-595 of human ER $\alpha$ .
<b>Clone:</b>	F3
<b>Isotype:</b>	IgG1, $\kappa$
<b>Application(s):</b>	ELISA, Western Blot, Immunoprecipitation, Immunocytochemistry (Formaldehyde fixation), Immunoprecipitation *, Ges shift *, Not Immunohistochemistry, ChIP-seq.  1/500-1/500 and *1/10-1/20
<b>Specificity:</b>	Does not cross react with chicken. Other reactivities not evaluated (see sequence homology)

### Product Overview

There are two classes of Estrogen Receptors (ER,) the first is intracellular and translocate to the Nucleus when activated through binding to 17 beta estradiol. They are named ER alpha and ER beta. The second G protein couples receptors which are membranes proteins and have intracellular signaling function. central ER $\alpha$  signaling is necessary for the actions of GLP-1 on food-reward behavior.

FP/18/20

#### For research laboratory use only - Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

#### Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac  
10 bis avenue Ampère  
78180 Montigny le Bretonneux - France  
Tel.: +33 (0)139 306 036  
<https://www.bertin-bioreagent.com/pa206/contact-us>  
<https://www.bertin-bioreagent.com>