



## RXR $\hat{\pm}$ (Retinoid X receptor $\hat{\pm}$ ) $\hat{A}$ Monoclonal Antibody

Cat No: G01034 - 100  $\hat{\mu}$ 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>	recombinant protein, epitope corresponding to DE-region of mouse RXR $\alpha$
<b>Clone:</b>	3A2
<b>Isotype:</b>	IgG1, $\kappa$



<b>Application(s):</b>	ELISA
	Western Blot
	Immunocytochemistry
	Immunohistochemistry
	Immunoprecipitation
	Gel shift
	Recommended dilutions: 1/500-1/5000

**Specificity:** mouse and human RXR $\alpha$ . No cross reactivity on RXR $\beta$  and RXR $\gamma$ .

### Scientific Literature

Rochette-Egly C, Lutz Y, Pfister V, Heyberger S, Scheuer I, Chambon P, Gaub MP. Detection of retinoid X receptors using specific monoclonal and polyclonal antibodies. Biochem Biophys Res Commun. 1994 Oct 28;204(2):525-36.

FP/20/20

For research laboratory use only  $\hat{c}$  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'Activit $\hat{e}$ s du Pas du Lac  
10 bis avenue Amp $\hat{e}$ re  
78180 Montigny le Bretonneux - France  
Tel.: +33 (0)139 306 036

<https://www.bertin-bioreagent.com/pa206/contact-us>  
<https://www.bertin-bioreagent.com>