



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

Cat No: G01035 - 100 $\hat{\mu}$ L

General Data

Shipping:	wet ice
Stability:	Store at -20degrees; shelf life 3 year maximum after production Avoid repeat freezing and thawing cycles. Does not contain any preservative
Formulation:	Liquid Ascite, does not contain any preservative therefore avoid repeat freezing and thawing cycles
Host:	Mouse
Antigen:	recombinant protein, epitope corresponding to DE-region of mouse RXR α .
Clone:	1D12
Isotype:	IgG1, κ
Application(s):	ELISA Gel shift (endogenous and recombinant protein)Recommended dilutions: 1/500-1/5000
Specificity:	mouse and human RXR α , RXR β and RXR γ .



Scientific Literature

Bortell R, Owen TA, Shalhoub V, Heinrichs A, Aronow MA, Rochette-Egly C, Lutz Y, Stein JL, Lian JB, Stein GS. Constitutive transcription of the osteocalcin gene in osteosarcoma cells is reflected by altered protein-DNA interactions at promoter regulatory

FP/01/20

For research laboratory use only $\hat{\text{c}}$ 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 $\hat{\text{e}}$ s du Pas du Lac
10 bis avenue Amp $\hat{\text{e}}$ 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>