



## CRABP II (Cellular retinoic acid binding protein II) Monoclonal Antibody

Cat No: G01004 - 100 µL

### 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
<b>Clone:</b>	4C9
<b>Isotype:</b>	IgG1, κ



<b>Application(s):</b>	ELISA Western Blot Immunocytochemistry recommended dilution: 1/500- 1/5000
------------------------	---

**Specificity:** mouse CRABP II. No cross reactivity on human CRABP II and on CRABP I.

### Product Overview

The vitamin A metabolite retinoic acid (RA) regulates gene transcription by activating several members of the nuclear receptor family of ligand-activated transcription factors: the classical RA receptors RAR $\alpha$ , RAR $\beta$ , and RAR $\gamma$  and the peroxisome proliferator-activated receptor  $\beta/\delta$  (PPAR $\beta/\delta$ ). The partitioning of the hormone between its receptors is regulated by two intracellular lipid-binding proteins, cellular retinoic acid-binding protein type II (CRABP-II), which delivers RA to RAR, and fatty acid-binding protein type 5 (FABP5), which shuttles it to PPAR $\beta/\delta$ .

### Scientific Literature

Gaub MP, Lutz Y, Ghyselinck NB, Scheuer I, Pfister V, Chambon P, Rochette-Egly C. Nuclear detection of cellular retinoic acid binding proteins I and II with new antibodies. J Histochem Cytochem. 1998 Oct;46(10):1103-11.

FP/19/20

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>