

Product Information

logo

CRABP I (Cellular retinoic acid binding protein I) Monoclonal Antibody

Cat No: G01002 - 100 µL

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:	synthetic peptide
Clone:	10F5
Isotype:	IgG1, κ
Application(s):	ELISA Western Blot Immunocytochemistry recommended dilution: 1/500-1/5000
Specificity:	mouse and human CRABPI. No cross reactivity on CRABPII.

Product Overview

Cellular retinoic acid-binding protein 1 is a protein that in humans is encoded by the CRABP1 gene. A number of specific carrier proteins for members of the vitamin A family have been discovered. CRABP is assumed to play an important role in retinoic acid-mediated differentiation and proliferation processes. CRABP1 is structurally similar to the cellular retinol-binding proteins, but binds only retinoic acid, and constitutively expressed. It is thought to play an important role in retinoic acid-mediated differentiation and proliferation processes. In addition, Crabp1 protects cells from excess RA by binding it in the cytosol, away from RARs.

Scientific Literature

Gaub MP, Lutz Y, Ghyselinck NB, Scheuer I, Pfister V, Chambon P, Rochette-Egly C. Nuclear detection of

FP/17/20

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>

Product Information

logo

CRABP I (Cellular retinoic acid binding protein I) Monoclonal Antibody

Cat No: G01002 - 100 µL

cellular retinoic acid binding proteins I and II with new antibodies. J Histochem Cytochem. 1998
Oct;46(10):1103-11.

FP/17/20

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>