

PRODUCT INFORMATION

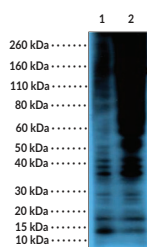
Phosphothreonine Monoclonal Antibody (Clone RM102)

Item No. 20716

Overview and Properties

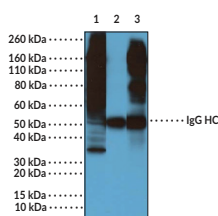
Contents:	This vial contains 100 µg of protein A affinity-purified antibody from an animal origin-free culture supernatant.
Immunogen:	Mixture of phosphothreonine-BSA conjugate and a phosphothreonine containing peptide
Cross Reactivity:	(+) Threonine-phosphorylated proteins, slight reactivity with phospho-serine peptides (-) Nonphosphorylated threonine, phosphoserine, phosphotyrosine
Species Reactivity:	(+) All species
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	50% glycerol/PBS with 1% BSA and 0.09% sodium azide
Clone:	RM102
Host:	Rabbit
Isotype:	IgG
Applications:	Chromatin IP (ChIP), ELISA, Flow cytometry (FC), Immunocytochemistry (ICC), Immunohistochemistry (IHC), Immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution is 1:100 to 1:500 for ChIP, ICC, IHC, and IP and 1:500 to 1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



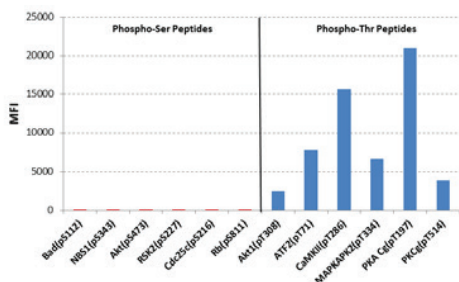
Lane 1: A431 (-)
Lane 2: CalA/Oka (+)

Western blot of serum-starved A431 cells nontreated or treated with Calyculin A/Okadaic Acid, using Phosphothreonine Monoclonal Antibody (Clone RM102) at a dilution of 1:2,000.

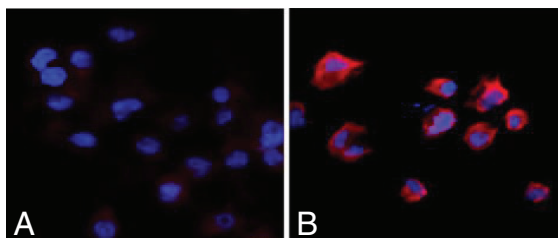


Lane 1: Whole lysate control
Lane 2: IP by rabbit IgG control
Lane 3: IP by RM102

Immunoprecipitation of Calyculin A/Okadaic Acid treated A431 cells by Phosphothreonine Monoclonal Antibody (Clone RM102) at a dilution of 1:500 and then blotted with RM102.



Phosphothreonine Monoclonal Antibody (Clone RM102) recognizes phosphorylated threonine in peptides with different sequences. It has minimal cross-reactivity with phosphorylated serine.



Panel A: Immunocytochemistry of serum-starved A431 cells nontreated or Panel B: Treated with Calyculin A/Okadaic Acid using Phosphothreonine Monoclonal Antibody (Clone RM102) at a dilution of 1:500 followed by a PE-conjugated secondary antibody, (red) and DAPI (blue).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/31/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Phosphothreonine Monoclonal Antibody (Clone RM102) is a probe for immunochemical detection of phosphorylated threonine residues on proteins by immunoblotting, immunohistochemistry, or immunocytochemistry. The phosphorylation of threonine residues by serine/threonine kinases serves a variety of purposes, including altering activity, stability, and interaction with other biomolecules. Phosphorylation may be persistent or transient, with dephosphorylation mediated by phosphatases.

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM