

# PRODUCT INFORMATION



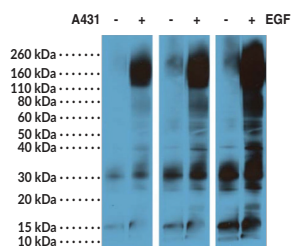
## Phosphotyrosine Monoclonal Antibody (Clone RM111)

Item No. 20717

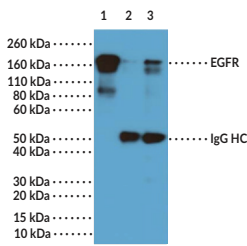
### Overview and Properties

<b>Contents:</b>	This vial contains 100 µg of protein A affinity-purified antibody from an animal origin free culture supernatant
<b>Immunogen:</b>	Phosphotyrosine-BSA conjugate
<b>Cross Reactivity:</b>	(+) Tyrosine-phosphorylated proteins; (-) Nonphosphorylated tyrosine, phosphoserine, phosphothreonine
<b>Species Reactivity:</b>	(+) All species
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	50% glycerol/PBS with 1% BSA and 0.09% sodium azide
<b>Clone:</b>	RM111
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	Western blot (WB), Chromatin IP (ChIP), ELISA, Flow Cytometry (FC), Immunocytochemistry (ICC), Immunohistochemistry (IHC), Immunoprecipitation (IP); the recommended starting dilution for WB is 1:1,000 to 1:5,000, 1:200 to 1:1,000 for IP and ChIP, and 1:200 to 1:500 for ICC and IHC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images

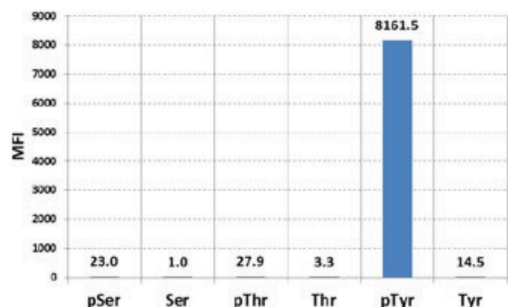


Western blot of serum-starved A431 cells nontreated or treated with EGF, using Phosphotyrosine Monoclonal Antibody (Clone RM111) at 1:5,000 dilution. The blot was exposed to the film from left to right at different time points.

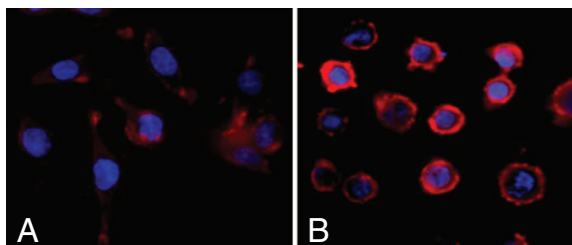


Lane 1: Whole lysate control  
Lane 2: IP by rabbit IgG control  
Lane 3: IP by Clone RM111

Immunoprecipitation of EGF-treated A431 cells by Phosphotyrosine Monoclonal Antibody (Clone RM111) 1:1,000 dilution, was blotted with an anti-EGFR rabbit monoclonal antibody.



Luminex analysis of Clone RM111's reaction against phosphorylated or nonphosphorylated serine, threonine, and tyrosine. Phosphotyrosine Monoclonal Antibody (Clone RM111) reacts only to phosphorylated tyrosine.



Panel A: Immunocytochemistry of serum-starved A431 cells nontreated or Panel B: Treated with EGF, using Phosphotyrosine Monoclonal Antibody (Clone RM111) at 1:500 dilution (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

---

Phosphotyrosine Monoclonal Antibody (Clone RM111) is a probe for immunochemical detection of phosphorylated tyrosine residues on proteins by immunoblotting. The phosphorylation of tyrosine residues by tyrosine 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