

Anti-RFP

Mouse Monoclonal Antibody

Catalog # R46-61M

Lot # W365-5

Cited Applications

- Western blot (1:1000)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes native and denatured forms of RFP protein.

Cross Reactivity

- Western blot RFP and its variants tag-RFP, DsRed, mCherry and mOrange, etc

Other RFP variants may also be detectable

Host/Isotype/Clone#

Mouse, IgG1, RF5R

Immunogen

RFP from the Discosoma sea anemone N-terminal peptide-KLH conjugates

Formulation

10mM PBS, pH7.2

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

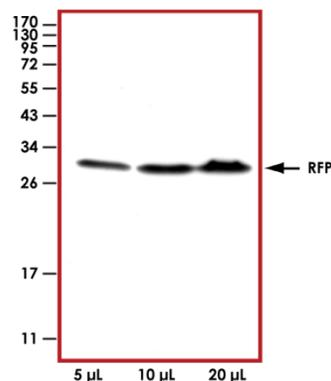
Scientific Background

RFP or red fluorescent protein from sea anemone *Entacmaea quadricolor* is a member of the green fluorescent protein family and is extensively used as a tag and reporter protein in molecular and cellular biology. RFP has a molecular weight of approximately 29 kDa and an excitation and emission maxima at 559 and 611 nm (1). RFP fluoresces over a wide pH range (4-10) with high quantum yield. Like other anthozoan fluorescent proteins, RFP forms tetramers at physiological concentrations.

References

1. Wiedenmann, J. et al: Red fluorescent protein eqFP611 and its genetically engineered dimeric variants. *J Biomed Opt.* 2005 Jan-Feb;10(1):14003..

Sample Data



Representative western blot with Anti-RFP (1:1000) using 5 µl, 10 µl and 20 µl of 293 cell lysate transfected with a RFP protein.

Anti-RFP

Mouse Monoclonal Antibody

Catalog Number

R46-61M

Specific Lot Number

W365-5

Purification

Affinity chromatography

Concentration

1.0 µg/µL

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.