



Glowing products for science

RFP Polyclonal Rabbit Antibody

This item is discontinued.

Product Description

This item is discontinued.

References

Download a list of [CF® dye references](#).

This datasheet was generated on June 3, 2026 at 03:01:25 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/rabbit-anti-rfp/>



Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Clonality	Polyclonal
Antibody type	Primary, Tag Antibody
Concentration	1 mg/mL
Host species	Rabbit
Antibody reactivity (target)	RFP
Cross adsorption	Human, Mouse, Rat
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide, Lyophilized: PBS/15 mg/mL BSA/20 mg/mL trehalose after reconstitution
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.