

# Product Information

## R-PE Conjugated Antibodies

Catalog No.	Product Description
20352	Goat Anti-Mouse IgG (H+L)
20353	Goat Anti-Rabbit IgG (H+L)
20354	Goat Anti-Rat IgG (H+L), highly cross-adsorbed Min X Hu, Bv, Hs and Rb
20355	Goat Anti-Human IgG (H+L), highly cross-adsorbed Min X Bv, Hs and Ms
20357	F(ab') <sub>2</sub> fragment of Goat anti-Mouse IgG (H+L)
20358	F(ab') <sub>2</sub> fragment Goat Anti-Rabbit IgG (H+L)

Bv: bovine; Ch: chicken; Gt: goat; GP: Guinea pig; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; Sh: sheep; SHm: Syrian hamster; Sw: swine; Rt: rat

### Unit Size

Whole IgG: 200 uL, 1 mL solution

F(ab')<sub>2</sub> fragment: 200 uL, 500 uL solution

**Concentration:** 0.5 mg/mL in pH ~7.4 PBS containing 2 mg/ml bovine serum albumin (IgG-free and protease-free) and 0.05% sodium azide.

**Form:** Red solution

### Spectral Property

$\lambda_{\text{abs}}/\lambda_{\text{em}}$  (in pH 7.4 PBS buffer): 496, 546, 555 nm/576 nm

### Storage and Handling

Store at 4°C, protected from light. Storage at -20°C is not recommended for R-PE conjugates. Product is stable for at least 6 months from date of receipt when stored as recommended.

### Product Description

The products are prepared by labeling IgG (H+L) with R-phycoerythrin. R-PE belongs to the family of phycobiliproteins, which are derived from cyanobacteria and eukaryotic algae. R-PE has absorption maxima at 496, 546, and 565 nm respectively and emission at 578 nm. Because of the high extinction coefficient (~1,960,000) and high fluorescence quantum yield (~0.82), R-PE is by far the brightest dye for the 565 nm wavelength region.

For most fluorescent secondary antibody applications, a concentration of 1 - 10 µg/mL is sufficient; however, optimal conditions should be determined empirically.

## Other R-PE bioconjugates

Cat. No.	Product Name	Unit Size
29045-100uL	Annexin V, R-PE Conjugate	20 assays
29045-500uL	Annexin V, R-PE Conjugate	100 assays
29044-200uL	Streptavidin, R-PE Conjugate	200 uL
29044-1mL	Streptavidin, R-PE Conjugate	1 mL

## Related Products

Cat. No.	Product Name	Unit Size
40061-T	RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO, Trial Size (15-20 tests)	25 uL
23001	EverBrite™ Mounting Medium	10 mL
23002	EverBrite™ Mounting Medium with DAPI	10 mL
23003	EverBrite™ Hardset Mounting Medium	10 mL
23004	EverBrite™ Hardset Mounting Medium with DAPI	10 mL
23005	CoverGrip™ Coverslip Sealant	15 mL
22005	Mini Super <sup>HT</sup> Pap Pen 2.5 mm tip, ~400 uses	1 pen
22006	Super <sup>HT</sup> Pap Pen 4 mm tip, ~800 uses	1 pen
22015	Fixation Buffer	100 mL
22016	Permeabilization Buffer	100 mL
22017	Permeabilization and Blocking Buffer	100 mL
22010	10% Fish Gelatin Blocking Buffer	100 mL
22011	Fish Gelatin Powder	2 x 50 g
22014	30% Bovine Serum Albumin Solution	100 mL
22002	Tween®-20	50 mL

Please visit [www.biotium.com](http://www.biotium.com) to view our full selection of secondary antibodies and other conjugates featuring bright and photostable fluorescent CF™ dyes, plus hundreds of other fluorescent reagents and assay kits for life science research.

For research use only.  
TWEEN is a registered trademark of Uniqema Americas LLC.