Rabbit IgG Isotype control polyclonal antibody (DyLight488™



Ordering Information

Order Online »

ENZ-ABS298-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable for at least 1 year after receipt when stored as recommended.

Handling Do not freeze. Protect from light.

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry

Application Notes Negative control in assays where Rabbit IgG (DyLight™

488)-labeled antibodies are used.

Emission Maximum 518 nm

Excitation Maximum 493 nm

Formulation Liquid. In PBS containing 0.09% sodium azide.

Host Rabbit

Isotype IgG

Source Purified from rabbit serum.

Species Reactivity Species independent

Technical Info / Product Notes DyLight™ 488 has high fluorescence intensity over a

broad pH range (pH 4-9) and is more photostable than many other dyes. The high water solubility of DyLight allows a high dye-to-protein ratio to be achieved without causing precipitation of the conjugates. Bright green fluorescence (λabs (max): 493nm, λem (max): 518nm,

εmax: 70,000).

Last modified: May 29, 2024



info-

eu@enzolifesciences.com