

Product Information

Andy Fluor™ 488 Goat Anti-Rabbit IgG (H+L) Antibody

Catalog Number	Unit Size
L110E	2 × 2.5 mL

Storage upon receipt:

- 4°C
- Protect from light

Ex/Em: 495/520 nm

Quick Start Guide



1. Apply 2 drops of antibody solution to fixed and permeabilized cells labeled with primary antibodies



2. Incubate 30 minutes at RT



3. Wash twice with PBS or HBSS



4. Image cells