

## Rabbit Anti-Fluorescein [FITC] Antibody, Biotinylated

Catalog Number: 46770

Unit Size: 100 ug

### Product Details

---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

Expiration Date	12 months upon receiving
-----------------	--------------------------

### Chemical Properties

---

Appearance	Liquid
------------	--------

Molecular Weight	N/A
------------------	-----

Soluble In	PBS
------------	-----

### Spectral Properties

---

Excitation Wavelength	N/A
-----------------------	-----

Emission Wavelength	N/A
---------------------	-----

### Applications

---

Rabbit Anti-Fluorescein [FITC] Antibody, Biotinylated is a polyclonal antibody that specifically binds to fluorescein-based dyes such as FITC and is directly labeled with biotin. This configuration enables flexible detection using streptavidin- or avidin-based systems in a variety of assay formats.

The antibody may cross-react with structurally related dyes such as Alexa Fluor® 488, DyLight® 488, and other fluorescein analogs. It is ideal for signal amplification in detecting low-abundance FITC conjugates, validating conjugation efficiency, and supporting multiplexed detection workflows.

*Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*