

## FITC (Fluorescein-5-isothiocyanate)

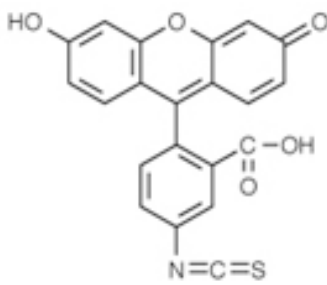
Catalog Number	Packaging Size
C100	50 mg
C101	500 mg

**Storage upon receipt:** -20°C, protect from light

### Introduction

**FITC** is a amine-reactive fluorescent probe for labeling of proteins. This amine-reactive fluorescein derivative has been the most common fluorescent derivatization reagent for covalently labeling of proteins.

### Specifications

<b>Label:</b>	Fluorescein	
<b>Ex/Em:</b>	495/518 nm	
<b>Detection Method:</b>	Fluorescent	
<b>Solubility:</b>	DMSO, DMF	
<b>Molecular Formula:</b>	C <sub>21</sub> H <sub>11</sub> NO <sub>5</sub> S	
<b>Molecular Weight:</b>	389.4	
<b>CAS Number:</b>	3326-32-7	
<b>Storage Conditions:</b>	-20°C, protect from light	
<b>Shipping Condition:</b>	Room Temperature	

### Applications

Protein labeling

### References:

1. Conjugation of antibodies with fluorochrome: modifications to the standard methods  
Goding, J.W.  
J Immunol Meth (1976) 13:215-226
2. Fluorescein labeling of Fab while preserving single thiol  
Der-Balian, G.P., et al.  
Anal Biochem (1988) 173:59-63
3. Fluorescence detection in automated DNA sequence analysis  
Smith, L.M., et al.  
Nature (1986) 321:674-678