



311PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# OneQuant™ Fluorescein Isothiocyanate (FITC)

(Cat. # 786-080)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

Fluorescein Isothiocyanate (FITC) is the simplest and most commonly used fluorescent dye. The isothiocyanate group of FITC will cross-link with amino, sulfhydryl, imidazolyl, tyrosyl or carbonyl groups on a protein. However, only the derivatives of primary and secondary amines generally yield stable products.

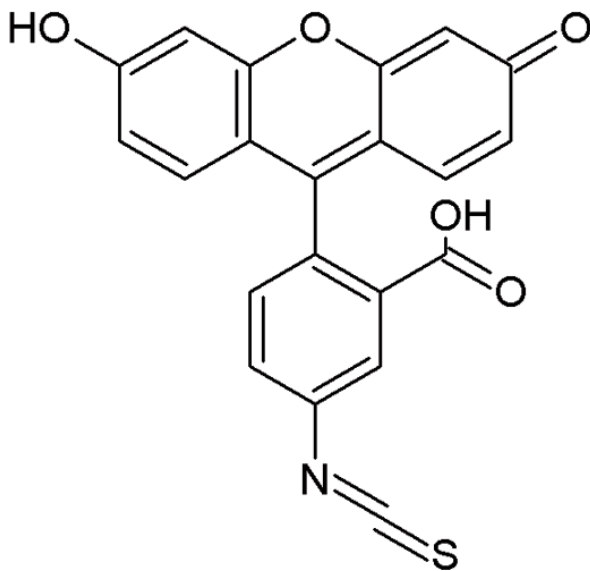
Chemical name: fluorescein-5-isothiocyanate

## SPECIFICATIONS

Formula:  $C_{21}H_{11}NO_5S$

Molecular Weight: 389.4

CAS No: 3326-32-7



## ITEM(S) SUPPLIED (Cat. # 786-080)

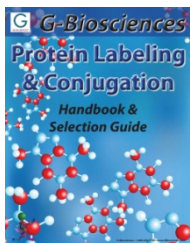
Description	Size
OneQuant™ Fluorecein Isothiocynate (FITC)	1mg/vial, 8 vials

## STORAGE CONDITIONS

It is shipped at ambient temperature. Upon receipt store product at  $-20^{\circ}\text{C}$  protected from moisture and light. Stable for one year.

## RELATED PRODUCTS

Download our Protein Labeling and Conjugation Handbook.



<http://info.gbiosciences.com/complete-protein-labeling-conjugation-handbook>

For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.

Last saved: 9/25/2012 CMH



[www.GBiosciences.com](http://www.GBiosciences.com)