

PRODUCT INFORMATION

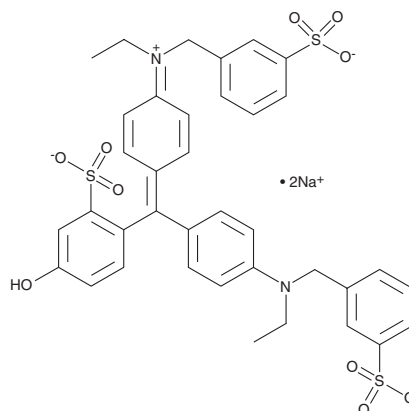


Fast Green FCF

Item No. 14335

CAS Registry No.: 2353-45-9
Formal Name: N-ethyl-N-[4-[[4-[ethyl[(3-sulfo-phenyl)methyl]amino]phenyl](4-hydroxy-2-sulfo-phenyl)methylene]-2,5-cyclohexadien-1-ylidene]-3-sulfo-benzenemethanaminium, disodium salt

Synonym: NSC 379443
MF: C₃₇H₃₄N₂O₁₀S₃ • 2Na
FW: 808.9
UV/Vis.: λ_{max}: 202, 306, 438, 616 nm
Supplied as: A crystalline solid
Storage: Room temperature
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Fast green FCF is supplied as a crystalline solid. A stock solution may be made by dissolving the fast green FCF in the solvent of choice, which should be purged with an inert gas. Fast green FCF is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of fast green FCF in ethanol and DMF is approximately 0.5 mg/ml and approximately 10 mg/ml in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of fast green FCF can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of fast green FCF in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Fast Green FCF is a dye used to stain proteins for IEF (isoelectric focusing) and SDS-PAGE. When bound to proteins, Fast Green FCF fluoresces near infrared (absorption maximum: 624 nm).¹ Fast Green FCF staining is linear over a wider range of protein concentrations compared to Brilliant Blue R.

Reference

1. Sabnis, R.W. Handbook of biological dyes and stains: Synthesis and industrial applications. John Wiley & Sons, Inc., Madison, NJ, USA (2010).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/14/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM