

# Product Information

## Methotrexate Conjugates

### Product List

Cat. No.	Conjugate	Ex/Em (nm)	MW
00091	CF@488A	490/516	1435
00092	CF@568	562/584	1234
00093	CF@594	593/615	1249
00094	CF@640R	642/663	1353
00023	Fluorescein, Triammonium Salt	498/517	979

**Unit Size:** 1 mg

### Storage and Handling

Store at -20°C. Protect from light. Product is stable for at least 12 months from date of receipt when stored as recommended.

Prepare the stock solution in DMSO and store at 4°C.

### Product Description

Methotrexate is a cytotoxic agent used as a chemotherapeutic and immunosuppressant. It is actively taken up by cells and binds to and inhibits dihydrofolate reductase, thereby inhibiting folic acid metabolism and inducing cell cycle arrest. Methotrexate fluorescein is useful for studying antimetabolite resistance and spontaneous gene amplification. Other dye conjugates of methotrexate may not be transported into cells, but may have other applications for studies of drug binding. Biotium offers methotrexate conjugated to fluorescein and a selection of our next-generation fluorescent CF@ Dyes.

Biotium's next-generation CF@ Dyes were designed to be highly water-soluble with advantages in brightness and photostability compared to other commercially available fluorescent dyes. For more information about CF@ Dyes, please visit our website at [www.biotium.com](http://www.biotium.com).

### References

1) J Cell Biol 162, 23 (2003); 2) Nat Biotechnol 19, 769-772 (2001); 3) J. Bol. Chem. 257, 14162(1982).

### Related Products

Cat. No.	Product
00060	Tetrodotoxin, Citrate Free
00061	Tetrodotoxin, With Citrate
00002... 00026	α-Bungarotoxin, CF@ Dye Conjugates
00068... 00079	Cholera Toxin Subunit B CF@ Dye Conjugates
00027... 00055	Phalloidin Conjugates
70030... 70034	Nerve Terminal Staining Kits
00081- 00087	Transferrin (Human) CF@ Dye Conjugates
22023	Paraformaldehyde, 4% in PBS Ready-to-Use Fixative
23001	EverBrite™ Mounting Medium
23002	EverBrite™ Mounting Medium with DAPI
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite™ Hardset Mounting Medium with DAPI
23008	Drop-n-Stain EverBrite™ Mounting Medium
23009	Drop-n-Stain EverBrite™ Mounting Medium with DAPI
23005	CoverGrip™ Coverslip Sealant
23023- 23024	Super HT Pap Pen 2.0
23012	TrueBlack® IF Background Suppressor System (Permeabilizing)
22010	10% Fish Gelatin Blocking Buffer
22014	30% Bovine Serum Albumin Solution
22002	Tween®-20
40046	Hoechst 33342, 10 mg/mL in H <sub>2</sub> O
40061	RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO

Please visit our website at [www.biotium.com](http://www.biotium.com) for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF@ Dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Tween is a registered trademark of Croda International LLC.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.