Transferrin (Human) CF® Dye Conjugates

Labeled human transferrin is used for microscopic studies of the endosomal pathway and/or transferrin uptake studies.

Diotium

Product Description

Labeled human holo-transferrin is used for microscopic studies of the endosomal pathway and/or transferrin uptake. Biotium's holo-transferrin conjugates are labeled with a selection our CFTM dyes, a series of next-generation fluorescent dyes developed at Biotium to have combined advantages in brightness, photostability, and water solubility compared to other fluorescent dyes.

- Choice of 7 CF® dyes from green to near-infrared
- Superior CF® dyes are bright, photostable, and water-soluble
- For endocytosis and transferrin uptake studies

Transferrin is a glycoprotein with homologous N-terminal and C-terminal iron-binding domains that binds up to two Fe(III) atoms for delivery to vertebrate cells through receptor-mediated endocytosis. After binding to its receptor on the cell surface, transferrin is rapidly internalized by invagination of clathrin-coated pits with formation of endocytic vesicles, the acidic environment favors dissociation of iron from the transferrin-receptor complex. Following the release of iron, the transferrin is then recycled back to the surface of the cell via the endocytic recycling pathway and released to bind more iron.

Superior CF® Dyes

Biotium's next-generation CF® dyes were designed to be highly water-soluble with advantages in brightness and photostability compared to Alexa Fluor®, DyLight®, and other fluorescent dyes. Learn more about CF® Dyes.

CF® Dye Human Transferrin

Product	Ex/Em	Size	Catalog No.
CF®488A Human Transferrin	490/515 nm	1 mg	Purchase 00081
CF®543 Human Transferrin	541/560 nm	1 mg	Purchase 00082
CF®568 Human Transferrin	562/583 nm	1 mg	Purchase 00083
CF®594 Human Transferrin	593/614 nm	1 mg	Purchase 00084
CF®640R Human Transferrin	642/662 nm	1 mg	Purchase 00085
CF®680R Human Transferrin	663/682 nm	1 mg	Purchase 00086
CF®750 Human	755/777 nm	1 mg	Purchase 00087

CF is a registered trademark of Biotium, Inc. Alexa Fluor, Texas Red, and DyLight are registered trademarks of Thermo Fisher Scientific.

References

Download a list of CF® dye references.

This datasheet was generated on November 5, 2025 at 12:03:33 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/human-transferrin-conjugate-cf-dye-conjugate/

Call us: 800-304-5357 Email: btinfo@biotium.com

Product attributes

Probe cellular localization	Membrane/vesicular	
For live or fixed cells	For live/intact cells	
Assay type/options	Real-time imaging	
Detection method/readout	Fluorescence microscopy, Flow cytometry	
Cell permeability	Membrane impermeant	
Fixation options	Fix before staining (formaldehyde), Fix after staining (formaldehyde), Permeabilize after staining	
Colors	Green, Orange, Red, Far-red, Near-infrared	