



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## ViaTag™ Haloalkane Ligands

Bright fluorescent ligands for the HaloTag® self-labeling protein tag. Available in 11 colors from blue to near-infrared, with options for labeling intracellular or cell surface targets.



### Product Description

ViaTag™ Haloalkane Ligands are fluorescent ligands that offer stable, covalent labeling of the HaloTag® self-labeling protein tag in live cells. They are offered in 11 different fluorescent colors from UV to near-IR, with both cell permeant and impermeant options available.

### Features

- Rapid covalent labeling of HaloTag® self-labeling tags
- Stable and ready-to-use DMSO stock solutions
- Minimal wash steps required
- Cell permeant and impermeant ligands, label live or fixed cells
- Wide selection of fluorescent colors from UV to near-IR
- Cost-effective alternative to other haloalkane ligands

The HaloTag® is a 34 kDa protein tag derived from bacterial haloalkane dehalogenase, which can be genetically fused to proteins of interest for expression in cells or organisms. The HaloTag® catalyzes the formation of a covalent bond with a fluorescent ligand containing a reactive haloalkane group, producing a stably labeled target protein detectable by fluorescence microscopy, flow cytometry, or other fluorescence-based methods.

Biotium's ViaTag™ Haloalkane Ligands take advantage of this covalent labeling system to provide bright, reliable fluorescence with minimal background. Available in both cell-permeant and cell-impermeant options, ViaTag™ ligands allow labeling of intracellular or cell surface HaloTag® fusion proteins in live or fixed cells. Because the labeling is covalent, it is resistant to fixation, permeabilization, and protein extraction. In addition, the tag enzyme retains activity after paraformaldehyde fixation, and can be labeled with ligand after fixation and permeabilization of the expressing cells.

ViaTag™ Haloalkane Ligands are offered in a broad selection of fluorescent colors and supplied as stable, ready-to-use DMSO stock solutions. Labeling is rapid and efficient, producing bright signal without the extensive wash steps required by other HaloTag® ligands.

### Choose from 11 bright colors, with flexible options for intracellular or cell surface targets

Biotium also offers a wide selection of [cellular stains](#) for labeling [membranes](#), [nuclei](#), and other organelles for live or fixed cells. This includes our [NucSpot® Live Cell Nuclear Stains](#) and [CellBrite® Steady Membrane Staining Kits](#) for long-term imaging of nuclei and membranes in live cells, respectively. Learn more about our unique [cellular stains](#) or view our [cellular stains guides](#) to view stains available for specific workflows or organisms.

### Product attributes

<b>For live or fixed cells</b>	Covalent & fixable stains, For live/intact cells
<b>Fixation options</b>	Fix after staining (formaldehyde), Permeabilize after staining
<b>Assay type/options</b>	Endpoint assay, Real-time imaging
<b>Colors</b>	Blue, Green, Red, Far-red, Near-infrared
<b>Storage Conditions</b>	Store at -10 to -35 °C, Protect from light

# ViaTag™ Haloalkane Ligands

Dye	Catalog No.	Ex/Em	Size	Cell membrane permeability	Laser line	Channel for detection
<a href="#">ViaTag™ 355/450 Haloalkane Ligand, 1000X in DMSO</a>	<a href="#">10301</a>	349/440 nm	30 uL	Permeant	355 nm (UV)	DAPI
<a href="#">ViaTag™ 400/460 Haloalkane Ligand, 1000X in DMSO</a>	<a href="#">10302</a>	391/452 nm	30 uL	Permeant	405 nm	DAPI, Alexa Fluor® 405, Pacific Blue®
<a href="#">ViaTag™ 500/525 Haloalkane Ligand, 1000X in DMSO</a>	<a href="#">10303</a>	504/527 nm	30 uL	Permeant	488 nm	FITC
<a href="#">ViaTag™ 540/565 Haloalkane Ligand, 1000X in DMSO</a>	<a href="#">10304</a>	543/572 nm	30 uL	Permeant	555 nm, 561 nm	Rhodamine, dsRed, PE
<a href="#">ViaTag™ 650/680 Haloalkane Ligand, 1000X in DMSO</a>	<a href="#">10305</a>	653/674 nm	30 uL	Permeant	640 nm	Alexa Fluor® 647, APC
<a href="#">ViaTag™ 490/525 Haloalkane Ligand, 1000X in DMSO</a>	<a href="#">10306</a>	493/517 nm	30 uL	Impermeant	488 nm	FITC
<a href="#">ViaTag™ 550/575 Haloalkane Ligand, 1000X in DMSO</a>	<a href="#">10307</a>	551/577 nm	30 uL	Impermeant	555 nm, 561 nm	Rhodamine, dsRed, PE
<a href="#">ViaTag™ 595/615 Haloalkane Ligand, 1000X in DMSO</a>	<a href="#">10308</a>	593/615 nm	30 uL	Impermeant	561 nm, 594 nm	Texas Red®
<a href="#">ViaTag™ 640/660 Haloalkane Ligand, 1000X in DMSO</a>	<a href="#">10309</a>	642/663 nm	30 uL	Impermeant	640 nm	Alexa Fluor® 647, APC
<a href="#">ViaTag™ 660/680 Haloalkane Ligand, 1000X in DMSO</a>	<a href="#">10310</a>	662/682 nm	30 uL	Impermeant	640 nm	Alexa Fluor® 660
<a href="#">ViaTag™ 740/765 Haloalkane Ligand, 1000X in DMSO</a>	<a href="#">10311</a>	742/767 nm	30 uL	Impermeant	730 nm	Alexa Fluor® 750

HALOTAG is a registered trademark of the Promega Corporation; ALEXA FLUOR, TEXAS RED, and PACIFIC BLUE are registered trademarks of Thermo Fisher Scientific.

This datasheet was generated on December 29, 2025 at 04:05:36 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/viatag-haloalkane-ligands/>