



Glowing products for science

SYT1 Recombinant Monoclonal Rabbit Antibody (rHL1626) - Biotium Choice

SYT1 (clone rHL1626) is a recombinant rabbit monoclonal antibody that recognizes Synaptotagmin 1 (SYT1). The antibody belongs to the Biotium Choice list of select antibodies that have been validated in-house for optimal performance.

Product Description

SYT1 Recombinant Monoclonal Rabbit Antibody (rHL1626) recognizes the Synaptotagmin 1 (SYT1) expressed in synaptic vesicles. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for fluorescence microscopy. The antibody is available conjugated to a selection of CF® Dyes. They are supplied in PBS, 2 mg/mL rBSA, and 0.05% sodium azide.

- Available in 8 bright and photostable CF® Dyes, including [near-IR CF®740](#) and [CF®790](#)
- Suitable for immunohistochemistry, immunofluorescence, and western

Synaptotagmin 1 (SYT1) belongs to the synaptotagmin membrane-trafficking protein family comprising 17 members. These proteins are single-pass transmembrane proteins characterized by two calcium-binding domains within their intracellular segment. SYT1 is predominantly found in presynaptic boutons of axon terminals, where it associates with synaptic vesicles. It is critical in initiating vesicle fusion and facilitating neurotransmitter release in response to elevated calcium levels in the local environment.

Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies - Powered by CF® Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes, they are our top-recommended antibodies for immunofluorescence and other applications.

Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for immunofluorescence and other applications
- Conjugated to bright, photostable [CF® Dyes](#) for superior signal and clarity
- New antibody clones and dye conjugates continuously in development

[View our full catalog of Biotium Choice antibodies](#)



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

Product attributes

Antibody number	#P055
Antibody reactivity (target)	SYT1
Antibody type	Biotium Choice, Primary
Biotium Choice Antibody	Primary
Host species	Rabbit
Clonality	Monoclonal, Recombinant Monoclonal
Clone	rHL1626
Isotype	IgG
Molecular weight	65-68 kDa
Synonyms	Synaptotagmin 1, P65, SVP65, SYT, AW124717, BAGOS, G630098F17Rik, P65, SVP65, SYT, SYT1, Sytl
Human gene symbol	SYT1
Entrez gene ID	6857
SwissProt	P21579
Immunogen	Recombinant fragment of human Synaptotagmin 1
Verified antibody applications	IF (verified), IHC (frozen) (verified)
Species reactivity	Human, Mouse, Rat
Expected antibody applications	IHC (FFPE) (published for clone), WB (published for clone)
Antibody application notes	Immunofluorescence (HeLa): 1-5 ug/mL Optimal concentration to be determined by end-user.
Positive control	Brain, retina
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: PBS/2 mg/mL rBSA/0.05% azide
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant rabbit IgG produced in Chinese Hamster Ovary (CHO) cell line
Antibody research areas	Neuroscience
Cell/tissue expression	Neurons

SYT1 Recombinant Monoclonal Rabbit Antibody (rHL1626) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	1 mL	P055-488A-1ML	CF®488A Features
CF®568	562/584 nm	100 ug/mL	1 mL	P055-568-1ML	CF®568 Features
CF®594	593/615 nm	100 ug/mL	1 mL	P055-594-1ML	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	1 mL	P055-640R-1ML	CF®640R Features
CF®647	652/668 nm	100 ug/mL	1 mL	P055-647-1ML	CF®647 Features
CF®680R	680/701 nm	100 ug/mL	1 mL	P055-680R-1ML	CF®680R Features
CF®740	742/767 nm	100 ug/mL	1 mL	P055-740-1ML	CF®740 Features
CF®790	783/808 nm	100 ug/mL	1 mL	P055-790-1ML	CF®790 Features

This datasheet was generated on August 21, 2026 at 01:08:15 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/syt1-recombinant-monoclonal-rabbit-antibody-rhl1626-biotium-choice/>