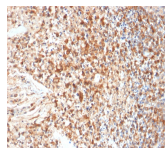


# Herpes Simplex Virus 1 (HSV1) Monoclonal Mouse Antibody (HSV1/2095)



## Product Description

The antibody reacts with HSV type 1 specific antigen. It is suitable for detection of HSV in human cellular material obtained from superficial lesions or biopsies and for the early identification of HSV in infected tissue cultures. The herpes simplex virus (HSV) (also known as cold sore, night fever or fever blister) is a virus that causes a contagious disease. There are two main types of Herpes Simplex Virus (HSV), 1 and 2. The HSV-1 strain generally appears in the orafacial organs. HSV2 usually resides in the sacral ganglion at the base of the spine. All herpes viruses are morphologically identical: they have a large double-stranded DNA genome and the virion consists of an icosahedral nucleocapsid, which is surrounded by a lipid bilayer envelope. UL42, the processivity subunit of the HSV-1 DNA polymerase binds DNA as a monomer and is essential for the replication of the virus. UL42 reduces the rate of dissociation from primer-template DNA, but it does not reduce the rate of elongation. UL42 increases the ability of UL9 to load onto DNA, thus increasing its assembly into a functional complex that is capable of unwinding duplex DNA. Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. **Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email [order@biotium.com](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

## Product attributes

Antibody number	#2095
Antibody reactivity (target)	HSV1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	HSV1/2095
Isotype	IgG1, kappa
Molecular weight	61 kDa
Synonyms	HSV1; Herpes simplex virus 1
Human gene symbol	Not Applicable
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
Unigene	Not Applicable
Immunogen	Baculovirus-expressed HSV DNA polymerase (POL) and POL/UL42 complex
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	HSV1
Positive control	HSV1 infected cells. Tissue.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, LI-COR Bioscience.