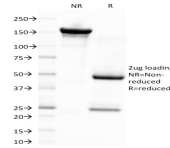


Epstein-Barr Virus Monoclonal Mouse Antibody (CS4)



Product Description

This antibody is a mixture of four different monoclonal antibodies. This antibody is specific to 60 kDa latent membrane protein (LMP-1) encoded by the BNLF1 gene of the EBV. Each clone reacts with different epitopes on the hydrophilic C-terminus of the cytoplasmic domain of LMP-1. This antibody stains strongly with EBV-positive lymphoblastoid cell lines and EBV infected B cell immunoblasts in infectious mononucleosis. EBV, also designated human herpesvirus 4 (HHV-4), is a member of the herpesvirus family and is one of the most common human viruses. EBV infects B cells and, though often asymptomatic, it can cause infectious mononucleosis, a disease characterized by fatigue, fever, sore throat and muscle soreness. 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 to inquire about stock status and lead times before placing your order.

Product attributes

Antibody number	#1744
Antibody reactivity (target)	Epstein-Barr Virus
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CS4
Isotype	IgG1, kappa
Molecular weight	60 kDa
Synonyms	EBNA1; EBV nuclear antigen 1; Epstein Barr nuclear antigen 1; Human Herpesvirus 4 (HHV4)
Human gene symbol	Not Applicable
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
Unigene	Not Applicable
Immunogen	Recombinant fusion protein containing the sequence of bacterial beta-galactosidase and the carboxyl half of EBV-encoded LMP
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	EBV
Positive control	EBV-infected cells. Tissues.
Shipping condition	Room temperature
Storage Conditions	Note: store BSA-free antibodies at -10 to -35°C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
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
BNC04	CF [®] 405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF [®] 488A	490/515	488	GFP, FITC
BNC68	CF [®] 568	562/583	532, 561	RFP, TRITC
BNC94	CF [®] 594	593/614	561	Texas Red [®]
BNC40	CF [®] 640R	642/662	633-640	Cy [®] 5
BNC47	CF [®] 647	650/665	633-640	Cy [®] 5
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

Dye Features

- [CF[®]405S Features](#)
- [CF[®]488A Features](#)
- [CF[®]568 Features](#)
- [CF[®]594 Features](#)
- [CF[®]640R Features](#)
- [CF[®]647 Features](#)

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.