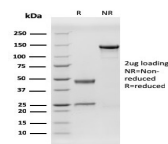


# Human Herpes Virus 8 Recombinant Monoclonal Mouse Antibody (HHV8/3606)



## Product Description

HHV 8 encodes a latent nuclear antigen (LNA), which is the product of the viral gene orf 73. LNA is capable of forming a complex with retinoblastoma susceptibility gene product, which may be related to its oncogenic activity. HHV8 is associated with three different diseases observed in AIDS patients; Kaposi's sarcoma, primary effusion lymphoma (which is a rare type of non-Hodgkin lymphoma affecting the body cavities) and multicentric Castlemann's disease. HHV 8 is the likely etiological agent of Kaposi sarcoma.

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.

**Catalog number key for antibody number 3606, Anti-Human Herpes Virus 8 (HHV8/3606)**

## Product attributes

<b>Antibody number</b>	#3606
<b>Antibody reactivity (target)</b>	Human Herpes Virus 8
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Recombinant Monoclonal
<b>Clone</b>	HHV8/3606
<b>Isotype</b>	IgG1, kappa
<b>Molecular weight</b>	Not Known
<b>Synonyms</b>	HHV8; Human Herpes Virus 8; Human herpesvirus 8; Kaposi's sarcoma associated herpes virus; KSHV
<b>Human gene symbol</b>	Not Applicable
<b>Entrez gene ID</b>	Not Applicable
<b>SwissProt</b>	Not Applicable
<b>Unigene</b>	Not Applicable
<b>Immunogen</b>	Recombinant protein corresponding to the latent nuclear antigen 1 molecule of HHV8
<b>Verified antibody applications</b>	IHC (FFPE) (verified)
<b>Species reactivity</b>	HHV8
<b>Antibody application notes</b>	ELISA: For coating use Ab at 1-5 ug/mL, order Ab without BSA; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody
<b>Positive control</b>	Herpes simplex type 1 (HSV-1) extract or infected cells. Tissue.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Product origin</b>	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
<b>Tumor expression</b>	Leukemia/lymphoma, Sarcoma

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>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 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, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on November 23, 2024 at 02:20:46 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/human-herpes-virus-8-recombinant-mouse-antibody-hhv8-3606/>