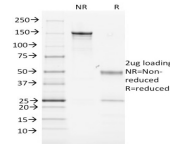


# MHC II, Mouse Monoclonal Mouse Antibody (MK-D6)



## Product Description

The antibody is directed against Mouse major histocompatibility complex class II, I-A d. It does not react with I-Ak, I-Ab, I-As, I-Af or I-Aa.

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.

## References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. PNAS USA (1982) 79:3604-3607. (functional studies)
2. PNAS USA (1987) 84: 2435-2439. (RIA)
3. J Inv Dermatol (1992) 99(5) 545-549. (Flow, surface; IF)
4. J Immunol (1996) 156: 4232-4239. (WB)

## Product attributes

<b>Antibody number</b>	#2007
<b>Antibody reactivity (target)</b>	MHC-II
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	MK-D6
<b>Isotype</b>	IgG2a, kappa
<b>Molecular weight</b>	kDa
<b>Synonyms</b>	Ia3; Ia-3; H-2Ea; H2-Ea; A1323765; I-Ealpha; E-alpha-f; MHC-H2-Ea
<b>Human gene symbol</b>	H2-Ea-ps (Mouse)
<b>Entrez gene ID</b>	100504404
<b>SwissProt</b>	P01904; P04224
<b>Unigene</b>	15680
<b>Immunogen</b>	Spleen cells from B10.D2 mice
<b>Antibody target cellular location</b>	Plasma membrane
<b>Species reactivity</b>	Mouse
<b>Expected antibody applications</b>	RIA (published for clone), Flow, surface (published for clone), Functional studies (published for clone), IF (published for clone), WB (published for clone)
<b>Positive control</b>	MPBMC's or Thymus.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
<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

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, and of LI-COR Bioscience.