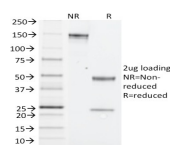


# NK1.1 / CD161c / Klr1c, Mouse Monoclonal Mouse Antibody (PK136)



## Product Description

NK1.1 is a type II integral membrane glycoprotein with a C-type lectin domain and is encoded by the *Klr1c/NKR-P1C* gene. It is predominantly expressed as a disulfide-linked homodimer on NK cells however; it is also expressed on NK-T cells, a rare population of T lymphocytes. NK1.1 is expressed in some mouse strains, including C57BL/6, FVB/N, and NZB, but not AKR, BALB/c, CBA/J, C3H, DBA/1, DBA/2, NOD, SJL, and 129. Expression of NKR-P1C antigen has been correlated with lysis of tumor cells in vitro and rejection of bone marrow allografts in vivo. NK-1.1 has also been shown to play a role in NK cell activation, IFN- $\gamma$  production, and cytotoxic granule release. NK-1.1 is commonly used as a mouse NK cell marker.

Primary antibodies are available purified, or with a selection of fluorescent CF<sup>®</sup> dyes and other labels. CF<sup>®</sup> dyes offer exceptional brightness and photostability. See the [CF<sup>®</sup> Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF<sup>®</sup>405S and CF<sup>®</sup>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<sup>®</sup> 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 2010, Anti-CD161c|Klr1c,|NK1.1 (PK136)**

## Product attributes

<b>Antibody number</b>	#2010
<b>Antibody reactivity (target)</b>	CD161c, Klr1c, NK1.1
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	PK136
<b>Isotype</b>	IgG2a, kappa
<b>Molecular weight of antigen</b>	kDa
<b>Synonyms</b>	CD161 antigen-like family member C; CD161c; Killer cell lectin-like receptor subfamily B member 1C (Klr1c); Klr1f1; Lymphocyte antigen 55c (Ly55c); Lymphocyte antigen 59; Natural killer cell surface protein P1-40; NK1.1; NKR-P1 40; NKR-P1.9; NKR-P1C; NKR-P1
<b>Human gene symbol</b>	Klr1c (Mouse)
<b>Entrez gene ID</b>	17059
<b>SwissProt</b>	P27814
<b>Unigene</b>	480764
<b>Immunogen</b>	CD Mouse spleen and bone marrow cells enriched for NK-1+ cells
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Species reactivity</b>	Mouse
<b>Expected antibody applications</b>	Flow, surface (published for clone), Functional studies (published for clone)
<b>Positive control</b>	MPBMC's or Thymus or Lymph Node.
<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
<b>Product origin</b>	Product may contain either bovine serum albumin (BSA) from bovine serum ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
<b>Cell/tissue expression</b>	NK cells

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF <sup>®</sup> 405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF<sup>®</sup>405S Features</a>
BNC88	CF <sup>®</sup> 488A	490/515	488	GFP, FITC	<a href="#">CF<sup>®</sup>488A Features</a>
BNC68	CF <sup>®</sup> 568	562/583	532, 561	RFP, TRITC	<a href="#">CF<sup>®</sup>568 Features</a>
BNC94	CF <sup>®</sup> 594	593/614	561	Texas Red <sup>®</sup>	<a href="#">CF<sup>®</sup>594 Features</a>
BNC40	CF <sup>®</sup> 640R	642/662	633-640	Cy <sup>®</sup> 5	<a href="#">CF<sup>®</sup>640R Features</a>
BNC47	CF <sup>®</sup> 647	650/665	633-640	Cy <sup>®</sup> 5	<a href="#">CF<sup>®</sup>647 Features</a>
BNC74	CF <sup>®</sup> 740	742/767	633-685	775/50	<a href="#">CF<sup>®</sup>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.

## References

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

J Exp Med (1993) 178: 223-229. (functional studies; Flow, surface)

This datasheet was generated on June 17, 2026 at 01:52:30 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/nk1-1-cd161c-klr1c-mouse-monoclonal-mouse-antibody-pk136/>