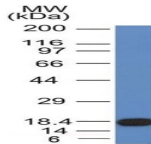


## CD117 / c-Kit Monoclonal Mouse Antibody (KIT/982)



### Product Description

CD117, also known as c-kit, is a member of the immunoglobulin superfamily. It is a receptor tyrosine kinase that is expressed on hematopoietic progenitor cells. It binds to the c-kit ligand KITLG/SCF and has roles in cell survival and cell proliferation. Upon ligand binding, it activates a number of signaling pathways. It is found on a wide variety of tumor cells including follicular and papillary carcinoma of thyroid, adenocarcinomas from endometrium, lung, ovary, pancreas, and breast as well as malignant melanoma, endodermal sinus tumor, and small cell carcinoma.

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 0982, Anti-CD117 (KIT/982)**

### Product attributes

Antibody number	#0982
Antibody reactivity (target)	CD117
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	KIT/982
Isotype	IgG1, kappa
Molecular weight of antigen	145 kDa
Synonyms	CD117; c-Kit; p145; Mast/stem Cell Growth Factor Receptor
Human gene symbol	KIT
Entrez gene ID	3815
SwissProt	P10721
Unigene	479754
Immunogen	Recombinant human CD117 fragment around aa100-300
Antibody target cellular localization	Plasma membrane
Verified antibody applications	WB (verified)
Species reactivity	Human
Antibody application notes	Immunohistology frozen only 2-4 ug/mL, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Western blotting 0.5-1.0 ug/mL, Optimal dilution for a specific application should be determined by user
Positive control	Esophagus, Gastrointestinal Stromal Tumor (GIST) or testicular germ cell tumor. Melanocytes in the basal layer of the epidermis and mast cells in the dermis of normal skin.
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Validated in protein array	Monospecific
Antibody research areas	Cancer, Signal transduction
Product origin	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.
Tumor expression	Breast cancer, Lung cancer, Melanoma, Ovarian cancer, Pancreatic cancer, Thyroid cancer

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 June 26, 2026 at 07:56:15 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/monoclonal-anti-cd117-kit982/>