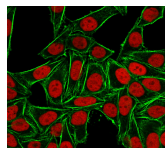


## Ku-Holo Monoclonal Mouse Antibody (KU729)



### Product Description

Recognizes a dimer of two proteins of 70 kDa and ~80 kDa, identified as two subunits of Ku. This MAB recognizes a conformational epitope of p70/p80 dimer, which is destroyed during Western blotting. The p70/p80 dimer is important for function of a 460 kDa DNA-dependent protein kinase. Ku protein plays a role in cell signaling, proliferation, DNA repair, replication, transcriptional activation, and apoptosis.

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 0729, Anti-Ku-Holo (KU729)**

### Product attributes

Antibody number	#0729
Antibody reactivity (target)	Ku-Holo
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	KU729
Isotype	IgG1, kappa
Molecular weight	70 kDa & 80 kDa
Synonyms	Ku (p70); 70kDa subunit of Ku antigen; ATP dependent DNA helicase 2 subunit 1; ATP-dependent DNA helicase II 70kDa subunit; CTC box-binding factor 75kDa subunit; CTC75; CTCBF; DNA repair protein XRCC6; G22P1; Ku autoantigen, 70kDa; Ku70; Kup70; Lupus Ku autoantigen protein p70; ML8; Thyroid autoantigen 70kD (Ku antigen); Thyroid-lupus autoantigen (TLAA); X-ray repair cross-complementing protein 6 (XRCC6) Ku (p80); 86kDa subunit of Ku antigen; ATP dependent DNA helicase 2 subunit 2; ATP dependent DNA helicase II 86kDa subunit; ATP-dependent DNA helicase II 80kDa subunit; CTC box-binding factor 85kDa subunit; CTC85; CTCBF; DNA repair protein XRCC5; KARP1; Ku autoantigen 80kDa; Ku80; Ku86 autoantigen related protein 1; KUB2; Lupus Ku autoantigen protein p86; Nuclear factor IV (NFIV); Thyroid-lupus autoantigen (TLAA); X-ray repair cross-complementing protein 5 (XRCC5)

Human gene symbol	XRCC6
Entrez gene ID	2547 (p70); 7520 (p80)
SwissProt	P12956 (p70); P13010 (p80)
Unigene	292493 (p70); 388739 (p80)
Immunogen	Nuclear extract of human HL-60 cells
Antibody target cellular	Nucleus
Verified antibody applications	Flow (intracellular) (verified), IF (verified)
Species reactivity	Human, Non-human primates
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunocytochemistry acetone-fixed cells 0.25-0.5 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Does not react with mouse, rat or chicken; others not known, Optimal dilution for a specific application should be determined by user
Positive control	Human cancer K562, HL-60, HEP-2 or HeLa cells
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
Antibody research areas	Apoptosis, Cell cycle, Signal transduction

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 Odyssey are registered trademarks of LI-COR Bioscience.

---

This datasheet was generated on August 29, 2024 at 08:27:41 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/monoclonal-anti-ku-holo-p7083-ku729/>