

# HL-60 Antigen / Human Leukemia Antigen Monoclonal Mouse Antibody (IPO-M6)

## Product Description

This antibody reacts with human leukemia cell line HL-60 and immunoprecipitates two proteins with MW of 48 kDa and 52 kDa. It does not stain B-cell lines Daudi, PHS, Namalwa, RPMI-1788 and T-cell lines CCRF-HSB2, Jurkat and Molt-4. This antibody stains monocytes and up to 10% of lymphocytes from peripheral blood of healthy donors. Blast cells of patients with AMMonL (M5 following FAB classification), AMMonL (M4) and hairy cells leukemia are positive. Its antigen is particularly expressed on blood cells from patients with infectious mononucleosis and CLL. Histiocytes and macrophages are also positive. Malignant cells from patients with AML (M1 and M2), T-ALL, B-ALL are not stained. HL-60 cells are used as an in vitro model of acute promyelocytic leukaemia and for differentiation and apoptosis studies. The HL60 cell line was established in 1977 from a patient with acute myeloid leukaemia. The cells largely resemble promyelocytes but can be induced to differentiate terminally in vitro. Some reagents cause HL60 cells to differentiate to granulocyte-like cells, others to monocyte/macrophage-like cells. The HL60 cell genome contains an amplified c-myc proto-oncogene; c-myc mRNA levels are correspondingly high in undifferentiated cells but decline rapidly following induction of differentiation. 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 0228, Anti-HL-60 Antigen (IPO-M6)**

## Product attributes

Antibody number	#0228
Antibody reactivity (target)	HL-60 Antigen, Human Leukemia Antigen
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	IPO-M6
Isotype	IgG1
Molecular weight	48 kDa & 52 kDa
Synonyms	Not Known
Human gene symbol	Not Known
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Leukemia HL-60 cells
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	Monocytes in tonsil or spleen
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	Immunology
Cell/tissue expression	Monocytes/macrophages
Tumor expression	Leukemia/lymphoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF®488A	490/515	488	GFP, FITC
BNC68	CF®568	562/583	532, 561	RFP, TRITC
BNC94	CF®594	593/614	561	Texas Red®
BNC40	CF®640R	642/662	633-640	Cy®5
BNC47	CF®647	650/665	633-640	Cy®5
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

## Dye Features

[CF®405S Features](#)
[CF®488A Features](#)
[CF®568 Features](#)
[CF®594 Features](#)
[CF®640R Features](#)
[CF®647 Features](#)

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 LI-COR Bioscience.