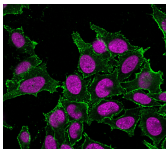




Glowing products for science

CD44 Standard Monoclonal Mouse Antibody (DF1485)



Product Description

Recognizes a cell surface glycoprotein of 80-95 kDa (CD44) on lymphocytes, monocytes, and granulocytes (Leucocyte Typing Workshop V). Its epitope is resistant to digestion by trypsin and chymotrypsin. The CD44 family of glycoproteins exists in a number of variant isoforms, the most common being the standard 85-95 kDa or hematopoietic variant (CD44s). Higher molecular weight isoforms are described in epithelial cells (CD44v), which are believed to function in intercellular adhesion and stromal binding. CD44 immunostaining is commonly used for the discrimination of urothelial transitional cell carcinoma in-situ from non-neoplastic changes in the urothelium.

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 to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1037, Anti-CD44 Standard (DF1485)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #1037 |
| Antibody reactivity (target) | CD44 Standard |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | DF1485 |
| Isotype | IgG1, kappa |
| Molecular weight | 80-95 kDa |
| Synonyms | LHR; BA-1; chondroitin sulfate proteoglycan 8 (CSPG8); Epican; Extracellular Matrix Receptor III (ECM III); GP90 Lymphocyte Homing Adhesion Receptor; HCAM; HCELL; hematopoietic cell E- and L-selectin ligand; Heparan Sulfate Proteoglycan; Hermes Antigen; HAS; HUTCH I; Hyaluronate Receptor; Indian blood group; Inlu Related p80 Glycoprotein; Ly 24; MDU2; MDU3; MIC4; MUTCH I; Phagocytic Glycoprotein 1 (PGP-1) |
| Human gene symbol | CD44 |
| Entrez gene ID | 960 |
| SwissProt | P16070 |
| Unigene | 502328 |
| Immunogen | Purified CD44 antigen (PGp-1) from lymphocyte membrane |
| Antibody target cellular localization | Plasma membrane |
| Verified antibody applications | IF (verified), IHC (FFPE) (verified) |
| Species reactivity | Human, Mouse |
| Expected antibody applications | Flow, surface (published for clone), IP (published for clone), WB (published for clone) |
| 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, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user |
| Positive control | HeLa cells or paracortex in tonsil or lymph node. |
| 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 |
| Product origin | Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |

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

Mol Cancer Ther (2011) 10(11): 2072-82. (western; immunoprecipitation; flow, surface)

This datasheet was generated on January 10, 2026 at 10:41:27 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/monoclonal-anti-cd44-standard-df1485/>

