

CD45 Mouse Monoclonal Antibody (2D1) - Biotium Choice

CD45 (clone 2D1) is a validated mouse monoclonal antibody that recognizes CD45, a transmembrane glycoprotein expressed in hematopoietic cells. The antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.



Product Description

CD45 (clone 2D1) is a mouse monoclonal antibody that recognizes transmembrane glycoprotein CD45. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance. The antibody is available conjugated to CF® Dyes and Astral™ tandem dyes. They are supplied in PBS, 0.1% BSA, 0.05% azide. CD45, also designated CD45R and PTPRC, has been identified as a transmembrane glycoprotein, broadly expressed among hematopoietic cells. Multiple isoforms of CD45R are distributed throughout the immune system according to cell type. These isoforms arise because of alternative splicing of exons 4, 5, and 6. The corresponding protein domains are characterized by the binding of monoclonal antibodies specific for CD45RA (exon 4), CD45RB (exon 5), CD45RC (exon 6) and CD45RO (exons 4 to 6 spliced out). The variation in these isoforms is localized to the extracellular domain of CD45R, while the intracellular domain is conserved. CD45R functions as a phosphor-tyrosine phosphatase. This MAb reacts with all isoforms of CD45R expressed by all hematopoietic cells, except erythrocytes, having a higher level of expression on lymphocytes than on granulocytes (Workshop IV). Antibody to CD45 is useful in differential diagnosis of lymphoid tumors from non-hematopoietic undifferentiated neoplasms. As an isotype control, we recommend [Isotype Control, Monoclonal Mouse IgG1 Kappa \(IGG1/1331\)](#) for this antibody. [View our full catalog of Biotium Choice antibodies](#), optimized and validated for the best results in flow cytometry.

Product attributes

Antibody number	P012
Antibody reactivity (target)	CD45
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Flow Antibodies, Primary
Host species	Mouse
Clonality	Monoclonal
Clone	2D1
Isotype	IgG1, kappa
Molecular weight	180 - 240 kDa (multiple isoforms)
Synonyms	Leukocyte Common Antigen (LCA), PTPRC, T200
Human gene symbol	CD45
Entrez gene ID	5788
SwissProt	P08575
Unigene	654514
Immunogen	Human
Antibody target cellular localization	Plasma membrane
Verified antibody	Flow (surface) (verified)
Species reactivity	Human
Antibody application notes	Recommended concentration for flow cytometry: 5 uL per million cells/0.1 mL
Positive control	Human peripheral blood mononuclear cells (PBMCs)
Shipping condition	Room temperature
Storage Conditions	Protect fluorescent conjugates from light
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Volume per assay	5 uL/test
Antibody/conjugate formulation	Conjugates: PBS/0.1% BSA/0.05% azide
Antibody research areas	Immunology
Cell/tissue expression	Lymphocytes

Conjugation	Ex/Em	Conc.	Size	Catalog No.
CF@405M	416/452 nm	20 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P012-405M-125 P012-405M-500
CF@488A	490/516 nm	200 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P012-488A-125 P012-488A-500
RPE-Astral™616	496, 546, 566/617 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P012-R616-125 P012-R616-500
CF@647Plus	652/668 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P012-647P-125 P012-647P-500
APC-Astral™813	633, 638/813 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P012-A813-125 P012-A813-500
R-PE	496, 546, 565/576 nm	50 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P012-RPE-125 P012-RPE-500
APC	651/660 nm	100 ug/mL	25 tests (125 uL) 100 tests (500 uL)	P012-APC-125 P012-APC-500

Dye Features

[CF@405M Features](#)
[CF@488A Features](#)