

## Human Lineage Marker/HLA-DR Antibody Mix (hLMAX)

**Catalog Number:** NBP2-29608

**Contents:** 25 tests of human FITC Lineage Markers & PerCP-Cy5.5 HLA-DR monoclonal antibody mix

**Formulation:** Tris or phosphate buffered saline with 0.05% sodium azide and 0.05% BSA

**Storage Conditions:** Store at 4°C, do not freeze; product is light-sensitive

**Test Size:** Novus' human Lineage Marker Antibody Mix contains 25 tests of each antibody, tested and standardized to use at 10 µl, to accommodate multi-color assay formats. Novus also offers service to undertake custom made conjugations with other fluorochromes, upon request.

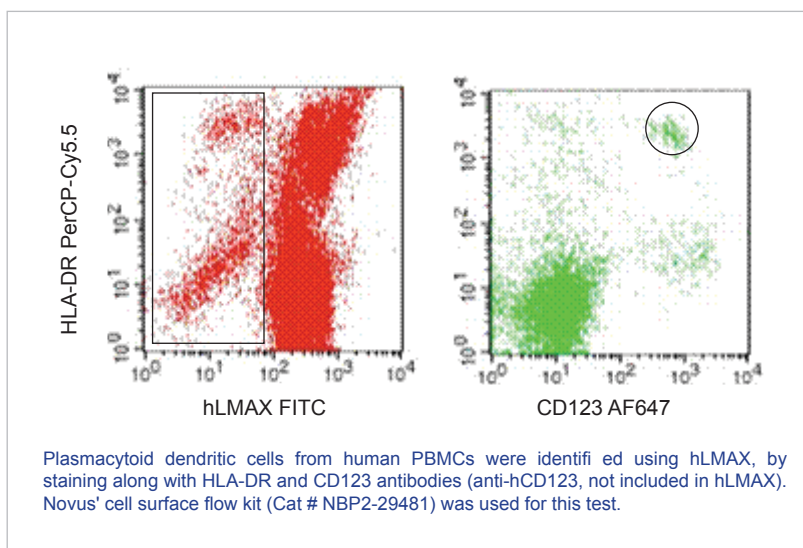
### Application Notes

Human LMAX (hLMAX) contains FITC-conjugated antibodies that, when added with HLA-DR as an additional marker, can be used to identify, enrich and/or deplete T, B, NK, and precursor myeloid cells (such as monocytes). The reagent can be used at 10 µl per one million cells in 100 µl total staining volume (or 100 µl whole blood). Novus' red blood cell lysis buffer (Cat # NBP2-29442) is recommended for lysing RBCs.

Human FITC Lineage Markers and PerCP-Cy5.5 HLA-DR monoclonal antibody mix has been validated in flow cytometric analysis of Plasmacytoid Dendritic Cells from:

- whole peripheral blood within 24h of collection
- fresh collected PBMCs
- frozen PBMCs

*This kit can also be used for the analysis of Myeloid Dendritic Cells.*



### hLMAX Components:

Antibody	Fluorochrome	Host	Isotype
CD3	FITC	Mouse	IgG2a
CD14	FITC	Mouse	IgG1
CD16	FITC	Mouse	IgG1
CD19	FITC	Mouse	IgG1
CD20	FITC	Mouse	IgG2b
CD56	FITC	Mouse	IgG1
HLA-DR	PerCP-Cy5.5	Mouse	IgG2a

### Product Citation:

#### **Tonsillar TLR9 expression and efficacy of tonsillectomy with steroid pulse therapy in IgA nephropathy patients.**

Sato D, Y Suzuki, T Kano, H Suzuki, J Matsuoka, H Yokoi, S Horikoshi, K Ikeda, Y Tomino. Nephrol Dial Transplant 0:1-8 (2011). **Novus product cited: Plasmacytoid Dendritic Cell (pDC)/TLR9 Kit (NBP2-29610). Flow (cell surface & intracellular): Primary human tonsillar B cells, Fig 3B. Note: NBP2-29608 is a component of the pDC/TLR9 (NBP2-29610) kit.**

---

For Research Use Only