

# PRODUCT INFORMATION



## HLA-DR Monoclonal Antibody (Clone L243)

Item No. 21827

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µg of protein G-purified IgG.
<b>Synonyms:</b>	HLA Class II Histocompatibility Antigen DR, MHC Class II Antigen DR
<b>Immunogen:</b>	Human lymphoblastoid cell line (RPMI 8866)
<b>Species Reactivity:</b>	(+) Human, canine, and non-human primate HLA-DR
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	PBS, pH 7.2, containing 50% glycerol, with 0.1% BSA, and 0.02% sodium azide
<b>Clone:</b>	L243
<b>Host:</b>	Mouse
<b>Isotype:</b>	IgG2a
<b>Applications:</b>	Flow cytometry (FC), Immunocytochemistry (ICC), Immunofluorescence (IF), Immunohistochemistry (IHC), Immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution for FC is 1:200. Suitable for ICC, IF, IHC, IP, and WB, working dilution should be determined empirically.

### Image

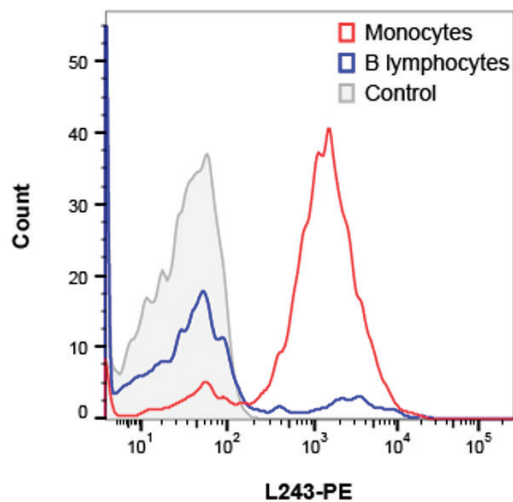


Figure 1: Flow cytometry data using PE conjugated HLA-DR Monoclonal Antibody (Clone L243). Highest levels detected in antigen presenting cell types.

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/21/2023

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

HLA-DR is a MHC Class II cell surface receptor heterodimer composed of a 33-35 kDa  $\alpha$  chain, a ~30 kDa  $\beta$  chain, and a 10-30 amino acid ligand.<sup>1</sup> When the heterodimer is fully combined on the cell surface of an antigen-presenting cell, such as macrophages, B cells, and dendritic cells, they present that ligand primarily to CD-4<sup>+</sup> T cells.<sup>2</sup> This presentation coupled with the T cell response can stimulate or suppress an antibody response to that ligand.<sup>2</sup> HLA-DRs have been linked to a number of autoimmune disorders such as rheumatoid arthritis, lupus, and psoriasis as well as diabetes, hepatitis, and sclerosis among a number of others.<sup>3-9</sup> Cayman's HLA-DR Monoclonal Antibody (Clone L243) detects HLA-DR with or without attached ligand. The predicated size of the HLA-DR $\alpha$  or HLA-DR $\beta$  subunits is ~29-30 kDa.

## References

---

1. Brodsky, F.M. and Parham, P. Monomorphic anti-HLA-A,B,C monoclonal antibodies detecting molecular subunits and combinatorial determinants. *J. Immunol.* **128(1)**, 129-135 (1982).
2. Denzin, L.K., Hammond, C., and Cresswell, P. HLA-DM interactions with intermediates in HLA-DR maturation and a role for HLA-DM in stabilizing empty HLA-DR molecules. *J. Exp. Med.* **184**, 2153-2165 (1996).
3. Panayi, G.S., Wooley, P., and Batchelor, J.R. Genetic basis of rheumatoid disease: HLA antigens, disease manifestations, and toxic reactions to drugs. *Br. Med. J.* **2(6148)**, 1326-1328 (1978).
4. Messner, R.P., De Horatius, R., and Ferrone, S. Lymphocytotoxic antibodies in systemic lupus erythematosus patients and their relatives: Reactivity with the HLA antigenic molecular complex. *Arthritis Rheum.* **23(3)**, 265-272 (1980).
5. Tsuji, K., Inouye, H., Nose, Y., et al. Further study on HLA-A, B, C, D, DR and haplotype antigen frequencies in psoriasis vulgaris. *Acta. Derm. Venereol. Suppl. (Stockh.)* **87**, 107-108 (1979).
6. Solow, H., Hidalgo, R., and Singal, D.P. Juvenile-onset diabetes HLA-A, -B, -C, and -DR alloantigens. *Diabetes* **28(1)**, 1-4 (1979).
7. Mackay, I.R. and Tait, B.D. HLA associations with autoimmune-type chronic active hepatitis: Identification of B8-DRw3 haplotype by family studies. *Gastroenterology* **79(1)**, 95-98 (1980).
8. Luckey, D., Bastakoty, D., and Mangalam, A.K. Role of HLA class II genes in susceptibility and resistance to multiple sclerosis: Studies using HLA transgenic mice. *J. Autoimmun.* **37(2)**, 122-128 (2011).
9. Gorodezky, C., Alaez, C., Murguía, A., et al. HLA and autoimmune diseases: Type 1 diabetes (T1D) as an example. *Autoimmun. Rev.* **5(3)**, 187-194 (2006).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM