

CD15 Recombinant monoclonal antibody (FUT4/1478R)

CD15 plays a role in mediating phagocytosis, bactericidal activity, and chemotaxis. It is present on >95% of granulocytes including neutrophils and eosinophils and to a lesser degree on monocytes. In addition, CD15 is expressed in Reed-Sternberg cells and some epithelial cells. CD15 antibody is very useful in the identification of Hodgkin's disease. CD15 is occasionally expressed in large cell lymphomas of both B and T phenotypes which otherwise have a quite distinct histological appearance.

This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

| | |
|-------------------------------------|---|
| Alternative Name | SSEA-1, Lewis X |
| Application | Flow Cytometry, IF, IHC (PS) |
| Clone | FUT4/1478R |
| Formulation | Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide. |
| Host | Rabbit |
| Immunogen | Purified human neutrophils. |
| Isotype | IgGk |
| Purity Detail | Protein A affinity purified. |
| Recommendation Dilutions/Conditions | Flow Cytometry (2-4 µg/106 cells in 0.1 mL)Immunofluorescence (2-4 µg/mL)Immunohistochemistry (formalin-fixed paraffin sections) (5-10 µg/mL for 30 min at RT)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. |
| Species Reactivity | Human, Mouse, Rat |
| Worry-free Guarantee | This antibody is covered by our Worry-Free Guarantee |