

PerCP Anti-human CD3 Antibody *OKT-3*Catalog number: 100341T0, 100341T1, 100341T2
Unit size: 25 tests, 100 tests, 500 tests**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.1 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse igg2a, κ
Immunogen	CD3e (T3E)
Clone	OKT-3
Conjugate	PerCP

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PerCP under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PerCP
Excitation Wavelength	477 nm
Emission Wavelength	678 nm

Applications

OKT-3 is an anti-human monoclonal antibody that targets the CD3e antigen. CD3e (alternatively called T cell antigen receptor complex or TCRC) is a 20 kD member of the Ig superfamily that is located on the surface of cells like T cells. CD3 is a member of important cellular pathways, in particular, the cell surface receptor signaling pathway, T cell receptor signaling pathway and negative regulation of smoothened signaling pathway. Furthermore, in certain organisms, it positively regulates calcium-mediated signaling, upregulates peptidyl-tyrosine phosphorylation and enhances cell-matrix adhesion. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands such as TCR. CD3 is a very popular antibody target, with over 80000 publications in the last decade. CD3e is vital to immunology research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was

purified through affinity chromatography and conjugated to PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 693/37 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant Analyzer 10).