

Anti-Dehydroepiandrosterone Antibody *Mouse anti-human, monoclonal IgG1* [V100205] [V100206]

Catalog Number:	V100205, V100206
Unit Size:	50 µg, 1 mg
Concentration:	1mg/ml
Molecular Weight:	150 Kd
Origin of Antibody:	Mouse
Other Names:	Dehydroepiandrosterone
Description:	Mouse anti-human, monoclonal IgG1
Purity:	>90% by SDS-Page and HPLC
Storage/Stability:	Store at -20°C/1 year. Avoid repeated freeze-thaw cycles
Form of Antibody:	Mouse IgG1 in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 15 mM sodium azide and 50% glycerol.
Applications:	ELISA