

Anti-Corticosterone antibody *Mouse anti-human, monoclonal IgG1* [V100210]

Catalog Number:	V100210, V100211	
Unit Size:	50 ug, 1mg	
Concentration:	1mg/ml	
Molecular Weight:	150 Kd	
Origin of Antibody:	Mouse ascites	
Other Names:	Corticosterone	
Description:	Mouse anti-human, monoclonal IgG1	
Purity:	>95% SDS-Page. Purified by ion exchange chromatography	
Specificity:	Corticosterone	100%
	17-beta-Estradiol	<3%
	Estrone	<5%
	Estriol	<6%
	17-alpha-Ethinylestradiol	<3%
	17-OH Progesterone	<1%
	Androstenedione	<1%
	Aldosterone	<1%
	17-OH Progesterone	<1%
	Progesterone	<1%
	Testosterone	<1%
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.	