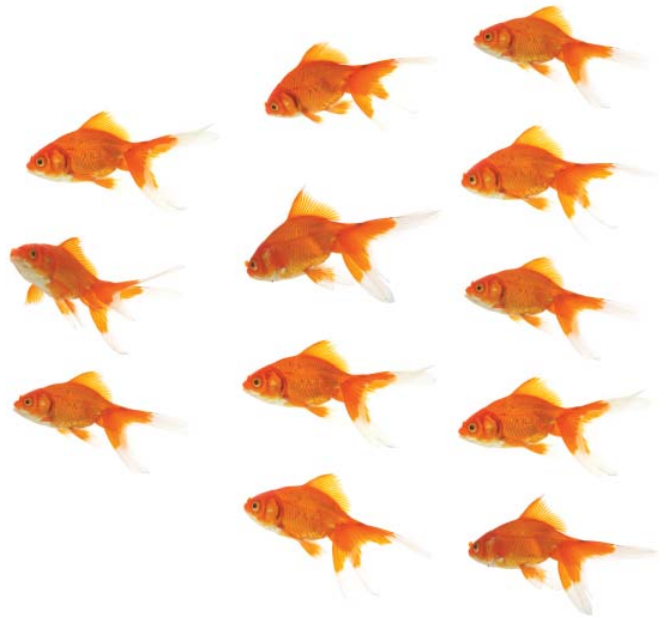


ELISA Kits for
MOG (1-125) & (35-55) IgG
Industry's First & Only



We're leading the way...

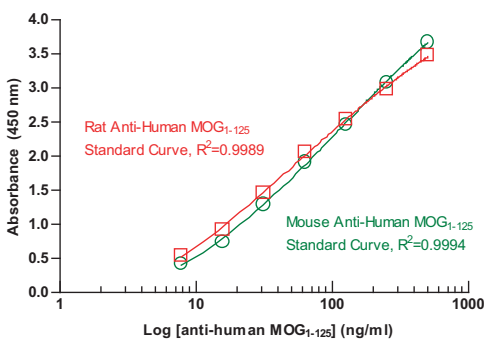


SensoLyte® Anti-MOG (1-125) & (35-55) IgG ELISA Quantitation Kits

AnaSpec's Anti-MOG ELISA assay kit series, the Sensolyte® Anti-Mouse MOG(1-125) IgG Quantitation Kit, Sensolyte® Anti-Rat MOG (1-125), Sensolyte® Anti-Human MOG (1-125) and Sensolyte® Anti-Mouse/Rat MOG (35-55) IgG ELISA assay kits are ideal for determining the amount of anti-mouse MOG (1-125), anti-rat MOG (1-125), anti-human MOG (1-125) and anti-mouse/rat MOG (35-55) antibodies, respectively. These kits, which represent the industry's first and only MOG ELISA kits, as well as the recombinant MOG (1-125) proteins are some of the examples of innovative products from AnaSpec.

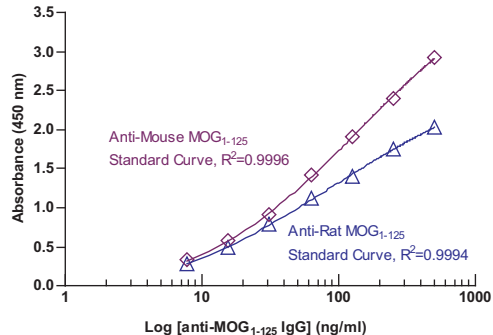
MOG (1-125) is found in the extracellular domain of Myelin Oligodendrocyte Glycoprotein (MOG). MOG, a member of the immunoglobulin superfamily, is expressed exclusively in the central nervous system (CNS). Although MOG protein constitutes only 0.01-0.05% of the CNS myelin proteins, it is implicated as a crucial autoantigen for multiple sclerosis in humans and experimental autoimmune encephalomyelitis (EAE) in rodents and monkeys.

4-PL Curve Fit for Rat and Mouse Anti-Human MOG₁₋₁₂₅ IgG Standards



Four parameter logistic (4-PL) curve fits for actual Mouse and Rat Anti-Human MOG₁₋₁₂₅ standards (kit component #55153B and #55153C).

4-PL Curve Fit for Anti-Rat and Anti-Mouse MOG₁₋₁₂₅ IgG Standards



Four parameter logistic (4-PL) curve fits for actual Anti-Mouse and Anti-Rat MOG₁₋₁₂₅ standards (kit components #55156B and #55157B).

MOG (1-125) & (35-55) IgG

ELISA Kits

Table 2. Comparison data of the Sensolyte® Anti-MOG IgG ELISA Quantitation Kits

	Sensolyte® Anti-Human MOG (1-125) Kit Cat# 55153	Sensolyte® Anti-Mouse MOG (1-125) Kit Cat# 55156	Sensolyte® Anti-Rat MOG (1-125) Kit Cat# 55157	Sensolyte® Anti-MOG (35-55) Kit (mouse/rat) Cat# 54465
Specificity	Human	Mouse	Rat	Mouse/Rat
Format	Human rMOG (1-125) precoated 12-strip plate	Mouse rMOG (1-125) precoated 12-strip plate	Rat rMOG (1-125) precoated 12-strip plate	MOG (35-55) precoated 12-strip plate
Standard(s) Included	Mouse & Rat anti-human MOG (1-125) IgG standards	Anti-mouse MOG (1-125) IgG standard	Anti-rat MOG (1-125) IgG standard	Anti-mouse/rat MOG (35-55) IgG standard
Sensitivity	Detects as low as 100 pg anti-human MOG (1-125)	Detects as low as 100 pg anti-mouse MOG (1-125)	Detects as low as 100 pg anti-rat MOG (1-125)	Detects as low as 100 pg anti-mouse/rat MOG (35-55)
Assay Time	2-3 hours assay time at room temperature			
Sample Concentration	0.5-1 µl of serum or cerebrospinal fluid			
Read-out	Colorimetric read-out at 450 nm absorbance			
Broad Dynamic Range	8-500 ng IgG/ml serum (depending on colorimetric developing time)			

Related Products

Table 3. Recombinant human, mouse and rat MOG proteins in 3 sizes.

Protein Species	Catalog #		
	100 µg	500 µg	1000 µg
Human	55158-100	55158-500	55158-1000
Mouse	55150-100	55150-500	55150-1000
Rat	55152-100	55152-500	55152-1000

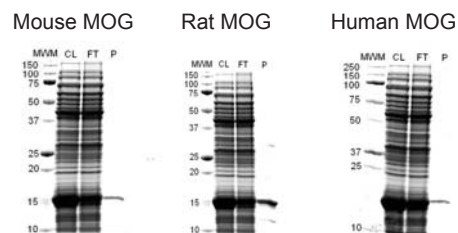


Fig. 1. purified recombinant MOG (rMOG) was loaded onto 10-20% Tris-HCl gel (10 µg/well) and resolved at 200V for 60 minutes. Protein markers and the purified rMOG (14.2 kDa) are indicated. CL= Crude Cell Lysate, FT= Flow Through, and P = Purified rMOG.

Table 4. Mono and polyclonal antibodies for MOG.

Antibody	Species Reactivity	Catalog#
Anti-MOG, mouse monoclonal clone B1A	Human, Mouse, Rat	55914
Anti-MOG, human, rabbit polyclonal	Human, Mouse, Rat	55161
Anti-MOG, mouse, rabbit polyclonal	Human, Mouse	55160

Table 5. Select listing of MOG peptides, with sequences from human (hu), mouse (ms) and/or rat (rt).

MOG (35-55), ms/rt Cat# 60130-1 (1 mg) Cat# 60130-5 (5 mg)	MOG (35-55), hu Cat# 63918 (1 mg)	MOG (37-54), ms/rt Cat# 62737 (1 mg)	MOG (38-55), ms/rt Cat# 61606 (1 mg)
MOG (43-54), hu, ms/rt Cat# 61610 (1 mg)	MOG (76-100), hu Cat# 64217 (1 mg)	MOG (91-114), rt Cat# 62738 (1 mg)	MOG (101-120), hu, ms Cat# 62370 (1 mg)



Certified to ISO 9001:2008