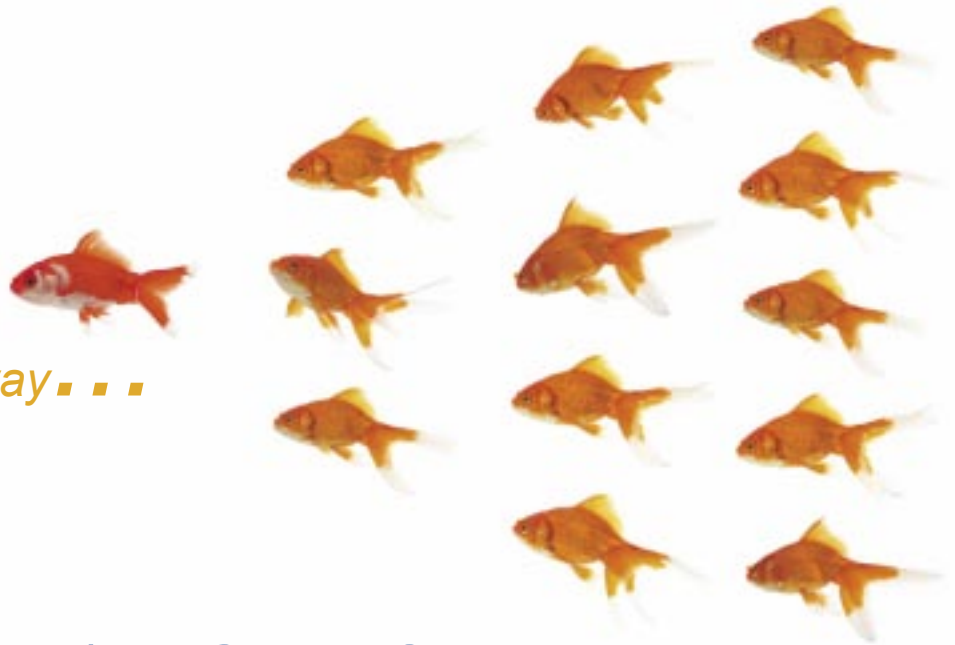


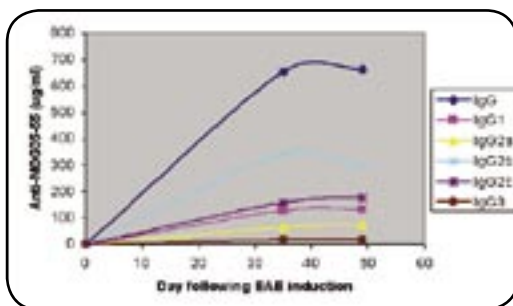
INNOVATION



We're leading the way. . . .

MOG (35-55) IgG ELISA Assay Kits Industry's First

Myelin oligodendrocyte glycoprotein (MOG) is a member of the immunoglobulin superfamily, which is expressed exclusively in central nervous system myelin. MOG (35-55) is able to induce autoantibody production and relapsing-remitting neurological disease with an extensive plaque like demyelination. The autoantibody response to MOG (35-55) has been seen in multiple sclerosis (MS) patients and MOG (35-55)-induced experimental autoimmune encephalomyelitis (EAE) mice. The exact pathological role of anti-MOG (35-55) autoantibody is under vigorous investigation. We have developed a series of useful and quantitative anti-MOG (35-55) ELISA kits for researchers to evaluate the role of anti-MOG (35-55) antibody in the development and treatment of EAE, an in vivo animal model for MS pathogenesis.



ISO 9001:2008 Company
AnaSpec, Inc. is a leading provider
of integrated proteomics solutions for
worldwide life science research.

MOG (35-55) IgG and Subtypes Quantitative Assay Kits

Convenient Format	Pre-coated and pre-blocked Ready to use (96-well strip plate) Ready to use substrate solution and other assay components
Minimal Sample	1-2 ul antiserum or body fluid
High Sensitivity	As little as 50 pg anti-MOG IgG
High Specificity	Extra-low cross reactivity when tested with the following rabbit antibodies Anti-PLP (139-151) @ 1:200 dilution Anti-PLP (178-191) @ 1:200 dilution Anti-MBP (1-11) @ 1:200 dilution
Broad Detection Range	0.5-200 ng antibody per ml serum (depending on colorimetric development time)

MOG (35-55) ELISA Kits	Catalog #
SensoLyte™ MOG (35-55) IgG Quantitative ELISA Kit (mouse, rat)	54465
SensoLyte™ MOG (35-55) IgG ₁ Quantitative ELISA Kit (mouse, rat)	54466
SensoLyte™ MOG (35-55) IgG _{2a} Quantitative ELISA Kit (mouse, rat)	54467
SensoLyte™ MOG (35-55) IgG _{2b} Quantitative ELISA Kit (mouse, rat)	54468
SensoLyte™ MOG (35-55) IgG _{2c} Quantitative ELISA Kit (mouse, rat)	54469
SensoLyte™ MOG (35-55) IgG ₃ Quantitative ELISA Kit (mouse, rat)	54470
SensoLyte™ MOG (35-55) IgG Subtype Quantitative ELISA Kit (mouse, rat)	54471

Related Products

MOG		MBP Peptides	
	Catalog #		Catalog #
Antibody			
Anti-MOG (35-55)	54003	MBP (1-11), rat: MBP (135-145), human	29200
Peptides		MBP (1-11), mouse	60627
MOG (35-55)	60130	MBP (1-17)	60625
MOG (38-53)	61613	[Ala4]-MBP (1-11)	29201
MOG (35-51)	61605	[Tyr4]-MBP (1-11)	29202
MOG (35-52)	61603	MBP (4-14)	22915
MOG (35-53)	61604	[Pyr4]-MBP (4-14)	23011
MOG (38-55)	61606	MBP (74-85), guinea pig	29207
MOG (40-54)	61626	[Ala81]-MBP (74-85)	29208
MOG (40-55)	61607	MBP (87-99), human	29211
MOG (41-54)	61608	[Arg91, Ala96]-MBP (87-99), human	29212
MOG (42-54)	61609	MBP (90-106)	60977
MOG (43-54)	61610	MBP (90-106), phosphorylated	60976
MOG (44-54)	61620	...more	
MOG (45-54)	61611		
MOG (46-54)	61612		
		PLP Peptides	Catalog #
		PLP (139-151)	29213
		[Ala144]-PLP (139-151)	29214
		[Gln144]-PLP (139-151)	29215
		[Leu144]-PLP (139-151)	29216
		[Leu144, Arg147]-PLP (139-151)	29217