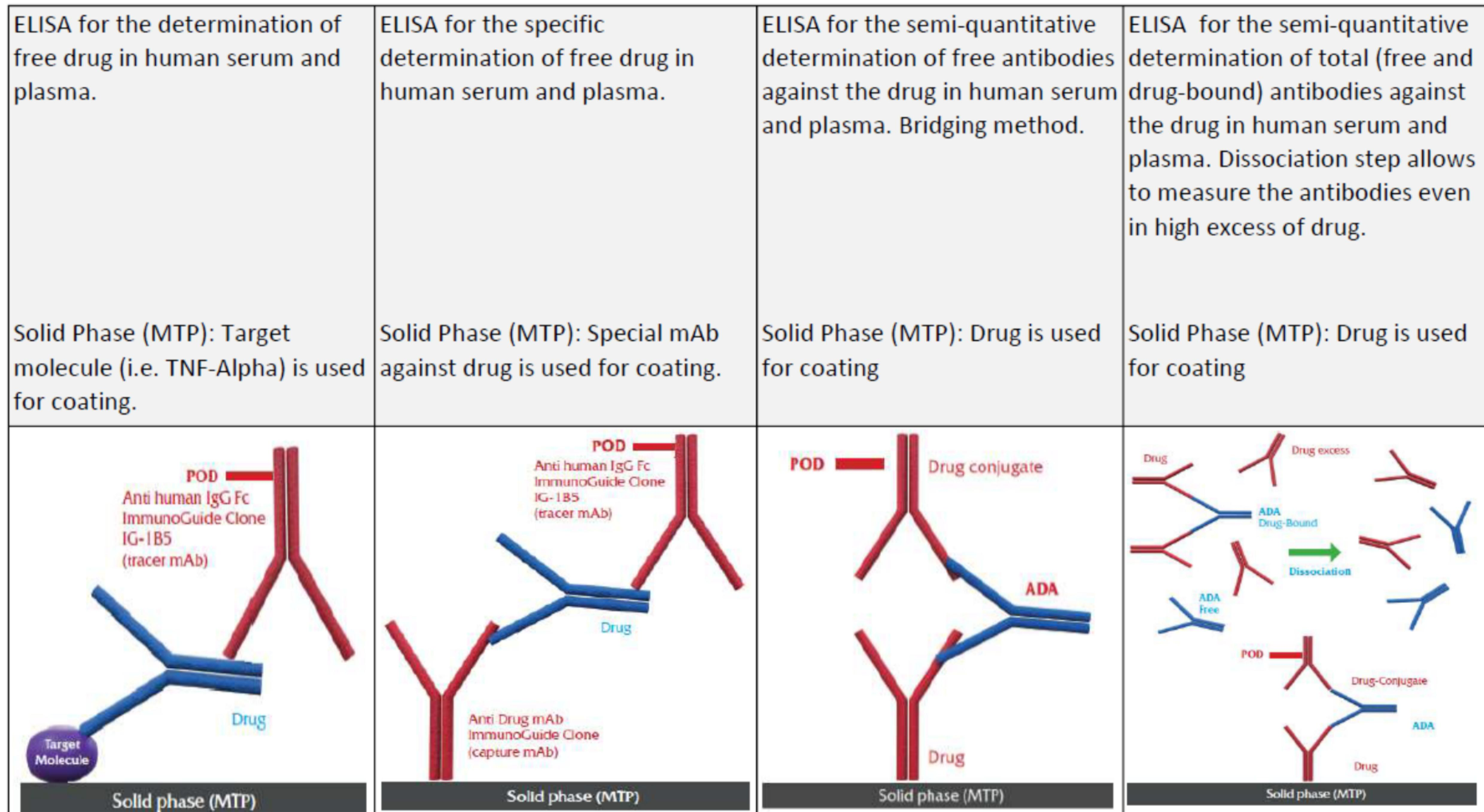




Drug Monitoring Assay Guide



Drug Name	Trade Name	Assay	mAb-Based Assay	Drug Antibodies Assay	Total Antibodies Assay
Infliximab	Remicade®, Remsima®, Inflectra®	IG-AA101	IG-AB101	IG-BB101	IG-TA101
Etanercept	Enbrel®	IG-AA102	IG-AB102	IG-BB102	
Adalimumab	Humira®	IG-AA103		IG-BB103	
Bevacizumab	Avastin®	IG-AA104	IG-AB104	IG-BB104	
Trastuzumab	Herceptin®, Herclon®	IG-AA105	IG-AB105	IG-BB105	
Rituximab	Rituxan®, Mabthera®		IG-AB106	IG-BB106	
Golimumab	Simponi®	IG-AA107	IG-AB107	IG-BB107	
Tocilizumab	Actemra®		IG-AB108	IG-BB108	
Certolizumab pegol	Cimzia®		IG-AB109	IG-BB109	
Omalizumab	Xolair®		IG-AB111	IG-BB111	
Cetuximab	Erbitux®		IG-AB112	IG-BB112	
Nivolumab	Opdivo®		IG-AB113	IG-BB113	
Aflibercept	Eylea®, Zaltrap®	IG-AA115			
Vendolizumab	Entyvio®		IG-AB116	IG-BB116	
Ustekinumab	Stelara®		IG-AB121		