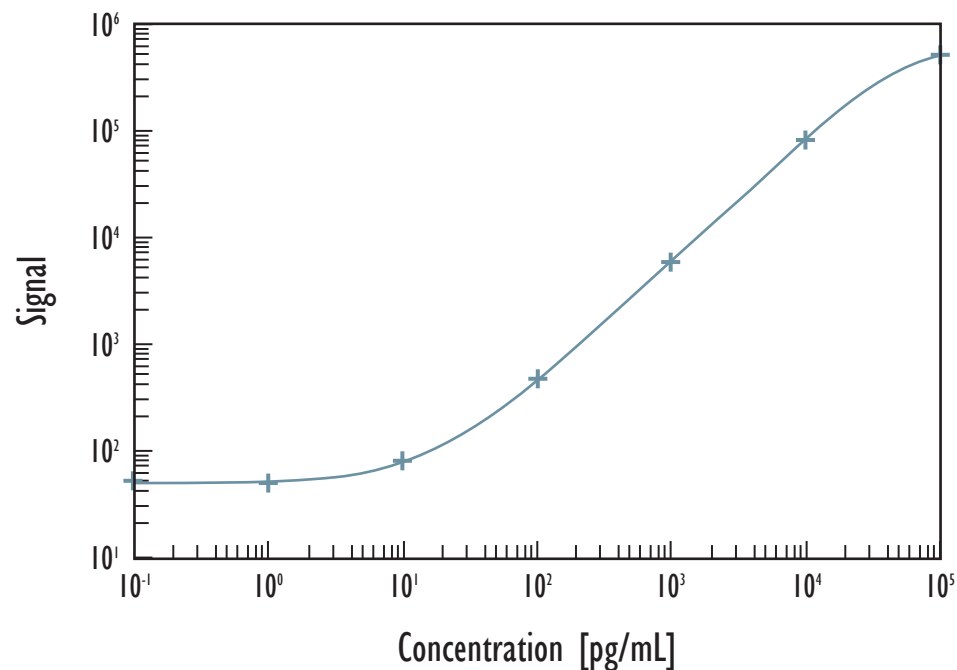


MULTI-ARRAY[®] HGF Assay

Detection of Hepatocyte Growth Factor in Serum and Plasma Samples



Sensitivity	10 pg/mL
Dynamic Range	10 to 100,000 pg/mL
Variability	Less than 10% CV

Spike Recovery and Dilution Linearity for Serum and Plasma Samples

	Spike or Dilution	Expected (pg/mL)	Measured (pg/mL)	% of Expected
Serum Pool	Neat		173	
	50% Dilution	87	86	100%
	Spike: +1,000 pg/mL	1173	1312	112%
EDTA Plasma Pool	Neat		119	
	50% Dilution	59	71	120%
	Spike: +1,000 pg/mL	1119	796	71%
Heparin Plasma Pool	Neat		76	
	50% Dilution	38	52	137%
	Spike: +1,000 pg/mL	1076	507	47%

Detection Limit	10 pg/mL
-----------------	----------

- Native HGF levels in pooled serum and plasma samples from apparently healthy donors are approximately tenfold above the detection limit of the HGF assay.
- The assay performs well for serum. Performance in EDTA plasma is better than in heparin plasma. (This may be because HGF contains a heparin-binding domain).

Kit Size	Catalog Number
1 plate	K151HDC-1
5 plates	K151HDC-2
20 plates	K151HDC-3
20 plates (Base)	K151HDA-3



Meso Scale Discovery
A division of Meso Scale Diagnostics, LLC.
www.mesoscale.com