



## Sample Reference Chart



= Validated

Kit/Assay/Analyte	Catalog #	Species	Serum	Plasma	TCM	Saliva	Tissue	Urine	Fecal	Citations
17-Hydroxyprogesterone	<a href="#">K053-H</a>	All								
20-Hydroxyecdysone	<a href="#">K066-H</a>	All								
Acetylcholinesterase (AChE) Activity Fluorescent	<a href="#">K015-F</a>	All								<a href="#">AChE Citations</a>
Aldosterone Chemiluminescent	<a href="#">K052-C</a>	All								
Aldosterone Colorimetric	<a href="#">K052-H</a>	All								<a href="#">Aldosterone Colorimetric Citations</a>
Allopregnanolone (monoclonal)	<a href="#">K061-H</a>	All								<a href="#">Allopregnanolone (monoclonal) Citations</a>
Allopregnanolone Chemiluminescent	<a href="#">K044-C</a>	All								
Allopregnanolone Colorimetric	<a href="#">K044-H</a>	All								<a href="#">Allopregnanolone Colorimetric Citations</a>
Arg8-Vasopressin Chemiluminescent	<a href="#">K049-C</a>	All								<a href="#">Arg8-Vasopressin Chemiluminescent Citations</a>
Atrial Natriuretic Peptide (ANP)	<a href="#">K026-H</a>	Mammal								
BCA Protein Colorimetric	<a href="#">K041-H</a>	All								<a href="#">BCA Protein Colorimetric Citations</a>
Butyrylcholinesterase (BChE) Fluorescent	<a href="#">K016-F</a>	All								<a href="#">BChE Fluorescent Citations</a>
Catalase Activity Colorimetric	<a href="#">K033-H</a>	All								<a href="#">Catalase Activity Colorimetric Citations</a>
Catalase Activity Fluorescent	<a href="#">K033-F</a>	All								<a href="#">Catalase Activity Fluorescent Citations</a>
Ceruloplasmin	<a href="#">K035-H</a>	All								<a href="#">Ceruloplasmin Citations</a>
Corticosterone Chemiluminescent	<a href="#">K014-C</a>	All								<a href="#">Corticosterone Chemiluminescent Citations</a>

Kit/Assay/Analyte	Catalog #	Species	Serum	Plasma	TCM	Saliva	Tissue	Urine	Fecal	Citations
Corticosterone Colorimetric	<a href="#">K014-H</a>	All								<a href="#">Corticosterone Colorimetric Citations</a>
Cortisol	<a href="#">K003-H</a>	All								<a href="#">Cortisol Citations</a>
Cortisone Chemiluminescent	<a href="#">K017-C</a>	All								<a href="#">Cortisone Chemiluminescent Citations</a>
Cortisone Colorimetric	<a href="#">K017-H</a>	All								<a href="#">Cortisone Colorimetric Citations</a>
Creatinine Serum	<a href="#">KB02-H</a>	Mammal								<a href="#">Creatinine Serum Citations</a>
Creatinine Serum Low Sample Volume	<a href="#">KB02-H1D</a>	Mammal								<a href="#">Creatinine Serum Low Sample Volume Citations</a>
Creatinine Urinary	<a href="#">K002-H</a>	All								<a href="#">Creatinine Urinary Citations</a>
Cyclic AMP Direct	<a href="#">K019-H</a>	All								<a href="#">Cyclic AMP Citations</a>
Cyclic AMP Direct Chemiluminescent	<a href="#">K019-C</a>	All								<a href="#">Cyclic AMP Direct Chemiluminescent Citations</a>
Cyclic GMP Direct	<a href="#">K020-H</a>	All								<a href="#">Cyclic GMP Direct Citations</a>
Cyclic GMP Direct Chemiluminescent	<a href="#">K020-C</a>	All								<a href="#">Cyclic GMP Direct Chemiluminescent Citations</a>
Cyclic GMP Direct Colorimetric	<a href="#">K065-H</a>	All								<a href="#">Cyclic GMP Direct Colorimetric Citations</a>
Cystatin C	<a href="#">K012-H</a>	Human								<a href="#">Cystatin C Citations</a>
Dehydroepiandrosterone Sulfate (DHEA-S)	<a href="#">K054-H</a>	All								
Direct 2'3'-Cyclic GAMP	<a href="#">K064-H</a>	All								
DNA Damage	<a href="#">K059-H</a>	All								
Endothelin-1 (ET-1)	<a href="#">K045-H</a>	Mammal								
Epiandrosterone	<a href="#">K063-H</a>	All								

Kit/Assay/Analyte	Catalog #	Species	Serum	Plasma	TCM	Saliva	Tissue	Urine	Fecal	Citations
Estrone	<a href="#">K031-H</a>	All								<a href="#">Estrone Citations</a>
Estrone-3-Glucuronide (E1G)	<a href="#">K036-H</a>	All								<a href="#">E1G Citations</a>
Estrone-3-Sulfate (E1S)	<a href="#">K038-H</a>	All								<a href="#">E1S Citations</a>
Formaldehyde Fluorescent	<a href="#">K001-F</a>	All								<a href="#">Formaldehyde Fluorescent Citations</a>
FRAP™ (Ferric Reducing Antioxidant Power)	<a href="#">K043-H</a>	All								<a href="#">FRAP Citations</a>
Galactose Colorimetric	<a href="#">K042-H</a>	All								
Glucose Colorimetric	<a href="#">K039-H</a>	All								<a href="#">Glucose Colorimetric Citations</a>
Glucose Fluorescent	<a href="#">K039-F</a>	All								
Glutathione Colorimetric	<a href="#">K006-H</a>	All								<a href="#">Glutathione Colorimetric Citations</a>
Glutathione Colorimetric Cuvette	<a href="#">K006-H1C</a>	All								<a href="#">Glutathione Colorimetric Cuvette Citations</a>
Glutathione Fluorescent	<a href="#">K006-F</a>	All								<a href="#">Glutathione Fluorescent Citations</a>
Glutathione Fluorescent 384-Well	<a href="#">K006-F1D</a>	All								<a href="#">Glutathione Fluorescent 384-Well Citations</a>
Glutathione Reductase Fluorescent	<a href="#">K009-F</a>	All								<a href="#">Glutathione Reductase Fluorescent Citations</a>
Glutathione S-Transferase Fluorescent	<a href="#">K008-F</a>	All								<a href="#">Glutathione S-Transferase Fluorescent Citations</a>
Histone Demethylase (HDM) Fluorescent	<a href="#">K010-F</a>	All								<a href="#">HDM Fluorescent Citations</a>
Hemoglobin Colorimetric	<a href="#">K013-H</a>	All								<a href="#">Hemoglobin Colorimetric Citations</a>
Hemoglobin High Sensitivity Colorimetric	<a href="#">K013-HX</a>	All								
Hydrogen Peroxide Colorimetric	<a href="#">K034-H</a>	All								<a href="#">Hydrogen Peroxide Colorimetric Citations</a>

Kit/Assay/Analyte	Catalog #	Species	Serum	Plasma	TCM	Saliva	Tissue	Urine	Fecal	Citations
Levonorgestrel (LNG)	<a href="#">K058-H</a>	All								
Myeloperoxidase Human	<a href="#">K060-H</a>	Human								
Nitric Oxide Colorimetric	<a href="#">K023-H</a>	All								<a href="#">Nitric Oxide Citations</a>
Osteopontin Human (OPN)	<a href="#">K021-H</a>	Human								
Oxytocin Chemiluminescent	<a href="#">K048-C</a>	All								
Oxytocin Colorimetric	<a href="#">K048-H</a>	All								<a href="#">Oxytocin Colorimetric Citations</a>
P450 Demethylation Activity	<a href="#">K011-F</a>	All								<a href="#">P450 Citations</a>
Palladium PdX™ API Fluorescent	<a href="#">K007-F</a>	All								
PGE2 Multi-Format Colorimetric	<a href="#">K051-H</a>	All								<a href="#">PGE2 Multi-Format Citations</a>
PGFM	<a href="#">K022-H</a>	All								<a href="#">PGFM Citations</a>
PKA (Protein Kinase A) Activity	<a href="#">K027-H</a>	All								<a href="#">PKA Citations</a>
Pregnanediol-3- Glucuronide (PDG)	<a href="#">K037-H</a>	All								<a href="#">PDG Citations</a>
Progesterone	<a href="#">K025-H</a>	All								<a href="#">Progesterone Citations</a>
Prolactin	<a href="#">K040-H</a>	All								
Retinol Binding Protein (RBP) (serum)	<a href="#">KU04-H</a>	All								<a href="#">RBP (serum) Citations</a>
Retinol Binding Protein (RBP) Multi-format	<a href="#">K062-H</a>	All								
Retinol Binding Protein High Sensitivity	<a href="#">KU04-H</a>	All								<a href="#">RBP High Sensitivity Citations</a>
ST2 Human	<a href="#">K055-H</a>									

Kit/Assay/Analyte	Catalog #	Species	Serum	Plasma	TCM	Saliva	Tissue	Urine	Fecal	Citations
Superoxide Dismutase (SOD) Colorimetric	<a href="#">K028-H</a>	All								<a href="#">Superoxide Dismutase Citations</a>
Testosterone	<a href="#">K032-H</a>	All								<a href="#">Testosterone Citations</a>
Thiol Fluorescent Detection	<a href="#">K005-F</a>	All								<a href="#">Thiol Fluorescent Detection Citations</a>
Thyroxine (T4)	<a href="#">K050-H</a>	All								<a href="#">T4 Publications</a>
Triiodothyronine (T3)	<a href="#">K056-H</a>	All								<a href="#">T3 Citations</a>
Urea Nitrogen (BUN)	<a href="#">K024-H</a>	All								<a href="#">Urea Nitrogen Citations</a>