

Steroid Hormone Assay Kits

Researchers have trusted Cayman's collection of steroid hormone assays for decades to study effects on metabolism, inflammation, immune function, reproduction, and stress responses.



- Assay Features:**
- Highly sensitive detection in the low pg/ml range
 - Validated in several biological matrices
 - Broad species reactivity

Androgens

Item No.	Product Name	Assay Range	Sensitivity (80% B/B ₀)
582701	Testosterone ELISA Kit	3.9-500 pg/ml	6 pg/ml
582751	11-keto Testosterone ELISA Kit	0.78-100 pg/ml	1.3 pg/ml
501740	Testosterone Glucuronide ELISA Kit	4.9-3,000 pg/ml	25 pg/ml

Estrogens

Item No.	Product Name	Assay Range	Sensitivity (80% B/B ₀)
501890	Estradiol ELISA Kit	0.61-10,000 pg/ml	20 pg/ml
582281	Estriol ELISA Kit	0.9-2,000 pg/ml	4 pg/ml (4°C for 18 hrs), 12 pg/ml (RT for 2 hrs)
501290	Estrone-3-Glucuronide ELISA Kit	2.7-6,000 pg/ml	23.2 pg/ml (overnight), 36.1 pg/ml (rapid)
582261	2-Methoxyestradiol ELISA Kit	13.1-8,000 pg/ml	40 pg/ml

Progestogens

Item No.	Product Name	Assay Range	Sensitivity (80% B/B ₀)
501300	Pregnanediol-3-Glucuronide (PDG) ELISA Kit	0.4-50 ng/ml	0.94 ng/ml
582601	Progesterone ELISA Kit	7.8-1,000 pg/ml	10 pg/ml

Glucocorticoids

Item No.	Product Name	Assay Range	Sensitivity (80% B/B ₀)
501320	Corticosterone ELISA Kit	8.2-5,000 pg/ml	30 pg/ml
500360	Cortisol ELISA Kit	6.6-4,000 pg/ml	35 pg/ml
500370	Cortisol Express ELISA Kit	39.1-5,000 pg/ml	110 pg/ml
501420	11-Oxoetiocholanolone ELISA Kit	0.061-1,000 ng/ml	0.2 ng/ml

Mineralocorticoids

Item No.	Product Name	Assay Range	Sensitivity (80% B/B ₀)
501090	Aldosterone ELISA Kit	15.6-2,000 pg/ml	30 pg/ml

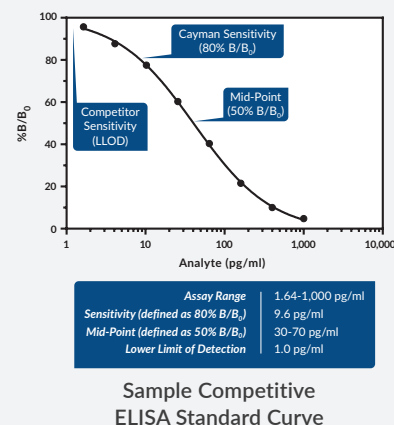
Vitamin D

Item No.	Product Name	Assay Range	Sensitivity (80% B/B ₀)
501050	Vitamin D ELISA Kit	0.19-25 ng/ml	0.5 ng/ml

Explore these products on our website for additional data, including mid-point (50% B/B₀) values and cross reactivity tables.

Why Choose Cayman's Steroid Assay Kits?

Cayman's steroid hormone immunoassays undergo rigorous quality testing to deliver the sensitivity and specificity needed to detect biologically significant analyte levels. These kits were developed to give the best sensitivity possible, while maintaining good signal-to-noise ratio, and report sensitivity as the 80% B/B₀ point on the standard curve. This point of quantitation can be trusted to give you reliable, reproducible data.



Find answers to frequently asked questions about Cayman's steroid hormone kits at www.caymanchem.com/steroidfaq

Steroid Hormone Receptors

Steroid hormones exert most of their effects by binding to and activating specific nuclear receptors, which are intracellular ligand-inducible transcription factors. Cayman offers cell-based assay kits to measure steroid hormone receptor activity in multiple species.

Androgen Receptor

Item No.	Product Name
15756	Human Androgen Receptor Reporter Assay System
28338	Rat Androgen Receptor Reporter Assay System
32578	Zebrafish Androgen Receptor Reporter Assay System

Estrogen Receptors (ERs)

ER α

Item No.	Product Name
15737	Human Estrogen Receptor Alpha Reporter Assay System
36709	Rat Estrogen Receptor Alpha Reporter Assay System
30256	Zebrafish Estrogen Receptor Alpha Reporter Assay System

ER β

Item No.	Product Name
15738	Human Estrogen Receptor Beta Reporter Assay System
36710	Rat Estrogen Receptor Beta Reporter Assay System

ER α & ER β

Item No.	Product Name
15739	Human Estrogen Receptors Reporter Assay Panel

Progesterone Receptor

Item No.	Product Name
15759	Human Progesterone Receptor Reporter Assay System
32581	Rat Progesterone Receptor Reporter Assay System

Glucocorticoid Receptor

Item No.	Product Name
15733	Human Glucocorticoid Receptor Reporter Assay System
21400	Mouse Glucocorticoid Receptor Reporter Assay System
33510	Rat Glucocorticoid Receptor Reporter Assay System
32579	Zebrafish Glucocorticoid Receptor Reporter Assay System

Mineralocorticoid Receptor

Item No.	Product Name
15740	Human Mineralocorticoid Receptor Reporter Assay System

Vitamin D Receptor

Item No.	Product Name
15742	Human Vitamin D Receptor Reporter Assay System



Manufactured by INDIGO Biosciences, Inc.

Explore all of our hormone and receptor assay kits
at www.caymanchem.com