

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

Adrenal Androgens & Androgen Precursor

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
502600	DHEA-S ELISA Kit	31.3-4,000 pg/ml	96 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
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
500370	Cortisol ELISA Kit (Express)	39.1-5,000 pg/ml	110 pg/ml
501420	11-Oxoetiocholanolone ELISA Kit	0.031-200 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



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

Steroid Hormone Receptors

Cell-based reporter assays complement ELISA by measuring functional receptor activation, confirming whether ligand binding translates into a biological response. They provide mechanistic insight into agonist/antagonist activity within a physiologically relevant cellular environment. 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

Progesterone Receptor

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

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

Thyroid Hormone Receptors (TRs)

TR- α

Item No.	Product Name
15749	Human Thyroid Hormone Receptor Alpha Reporter Assay System

TR- β

Item No.	Product Name
15750	Human Thyroid Hormone Receptor Beta Reporter Assay System
30258	Zebrafish Thyroid Hormone Receptor Beta Reporter Assay System

TR- α & TR- β

Item No.	Product Name
15751	Human Thyroid Hormone Receptor Panel

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