

VALIDATION DATA

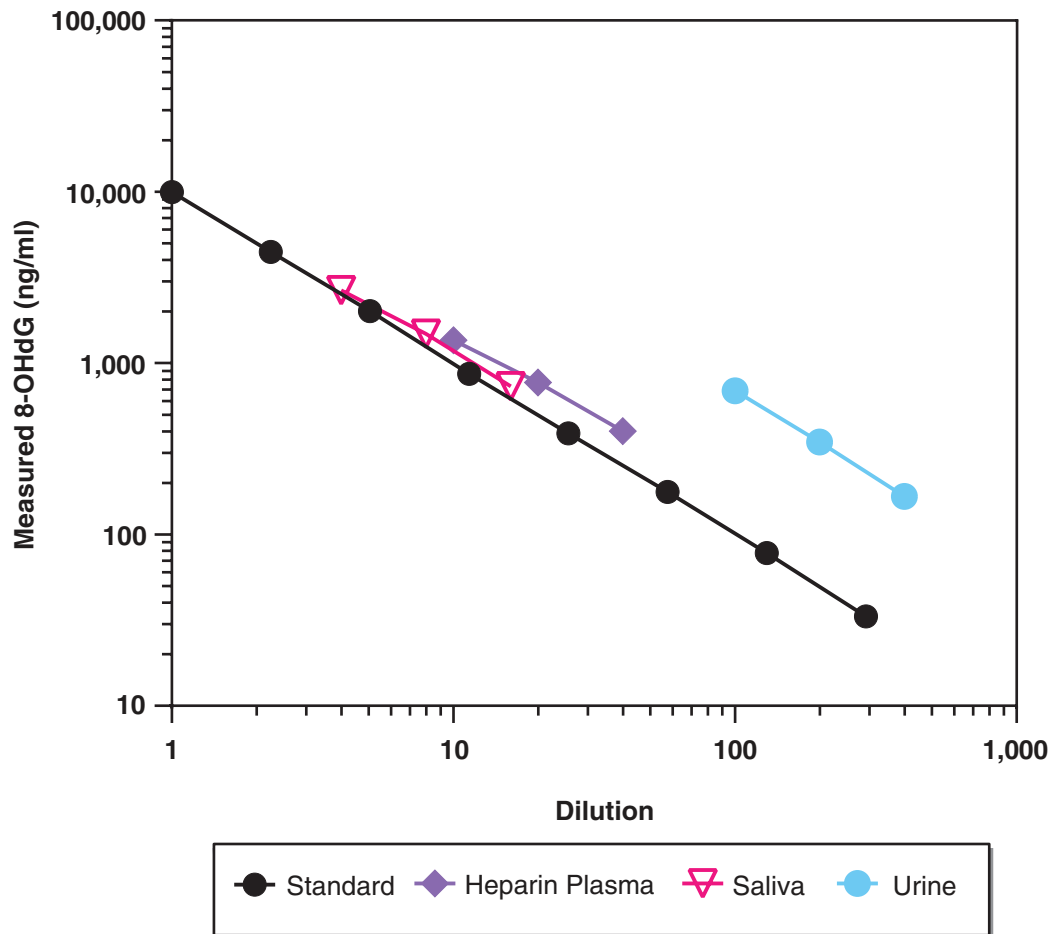


DNA/RNA Oxidative Damage Express ELISA Kit

Item No. 502890

Parallelism

To assess parallelism, human plasma, urine, and saliva and were assayed at multiple dilutions using the DNA/RNA Oxidative Damage Express ELISA. Concentrations were plotted as a function of sample dilution. The results are shown below.



WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/02/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

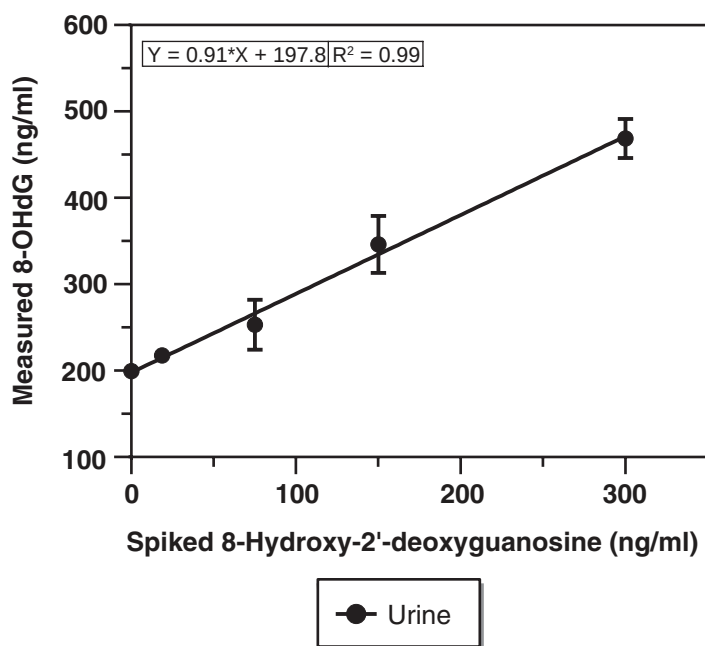
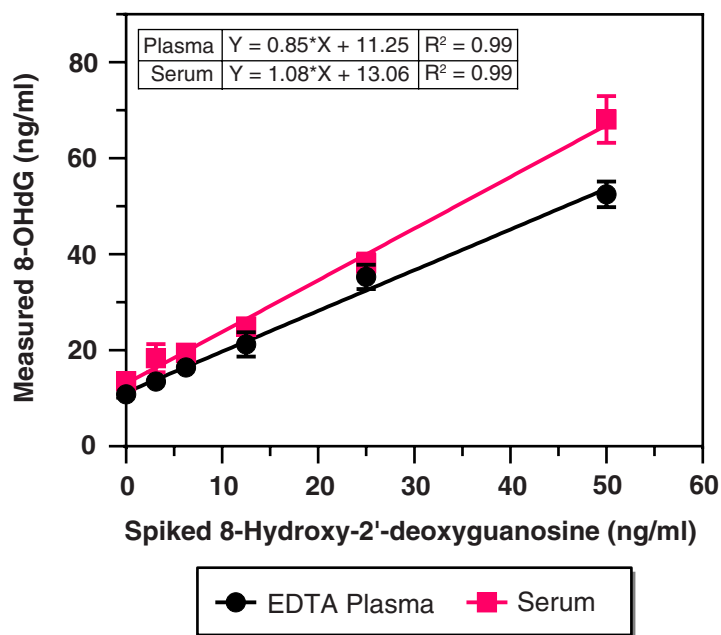
CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

VALIDATION DATA

Spike and Recovery

Human plasma, serum, and urine were spiked with 8-hydroxy-2'-deoxyguanosine (8-OHdG), processed as described in the Sample Preparation section (in the kit booklet), serially diluted with Assay Buffer, and evaluated using DNA/RNA Oxidative Damage Express ELISA. The results are shown below. The error bars represent standard deviations obtained from multiple dilutions of each sample.



CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

VALIDATION DATA



Linearity

Human plasma, serum, and urine samples were spiked with 8-OH-dG, processed as described in the Sample Preparation section, serially diluted with Assay Buffer, and evaluated for linearity using the DNA/RNA Oxidative Damage Express ELISA. The results are shown in the table below.

Dilution Factor	Measured Concentration (ng/ml)	Linearity%
EDTA Plasma, spiked with 25 ng/ml		
25	33.4	100
50	37.1	111
100	37.3	112
Serum, spiked with 25 ng/ml		
25	35.0	100
50	36.1	103
100	39.2	112
Urine, spiked with 300 ng/ml		
200	443	100
400	478	108
800	485	109

NOTE: Linearity has been calculated using the following formula:

$\% \text{Linearity} = (\text{Observed concentration value, dilution adjusted} / \text{First observed concentration value in the dilution series, dilution adjusted}) * 100$

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

VALIDATION DATA



Precision

Intra-assay precision was determined by analyzing 24 replicates of three controls in a single assay.

Matrix Control	Measured 8-OH-dG (ng/ml)	%CV
Control 1 (urine)	146	5.4
Control 2 (human heparin plasma)	51.3	6.9
Control 3 (human heparin plasma)	15.7	5.4

Inter-assay precision was determined by analyzing replicates of three controls in 8 separate assays on different days.

Matrix Control	Measured 8-OH-dG (ng/ml)	%CV
Control 1 (urine)	152	13.5
Control 2 (human heparin plasma)	50	5.9
Control 3 (human heparin plasma)	15.8	6.7

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM