

MR Bioassay for Environmental Sample Monitoring

The principal application of this Mineralocorticoid Receptor bioassay is to screen processed environmental samples, with the aim of quantifying the aggregate bioactivities, if any, of contaminating endocrine disrupting chemical (EDC) entities. This assay utilizes proprietary mammalian cells expressing the native human Mineralocorticoid Receptor (MR; nr3c2), a ligand-dependent transcription factor. The cells also include the luciferase reporter gene functionally linked to a MR-responsive promoter. Thus, quantifying changes in luciferase expression in the treated reporter cells provides a sensitive surrogate measure of the changes in total MR activity.

Ordering Information

[Order Online »](#)

NGO-EM00501-0096	96 wells
------------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping

Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name

Mineralocorticoid Receptor

Contents

This Mineralocorticoid Receptor Bioassay for Environmental Sample Monitoring kit is an all-inclusive luciferase reporter assay system that includes, in addition to MR Reporter Cells, two optimized media for use during cell culture and in diluting the user's test samples, the reference agonist Aldosterone, Luciferase Detection Reagent, a cell culture ready assay plate, and a detailed protocol and detailed protocol.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com