

# AR Bioassay for Environmental Sample Monitoring

The principal application of this assay is to functionally screen processed water samples, with the aim of quantifying the bioactivities, if any, of contaminating endocrine disrupting chemical (EDC) entities. This assay utilizes proprietary mammalian cells expressing the native Androgen Receptor (AR; NR3C4), a ligand-dependent transcription factor. The cells also include the luciferase reporter gene functionally linked to an AR-responsive promoter. Thus, quantifying changes in luciferase expression in the treated reporter cells provides a sensitive surrogate measure of the changes in total AR activity.

## Ordering Information

[Order Online »](#)

NGO-EM03001-0096	96 wells
------------------	----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Shipping

Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name**

Androgen Receptor

**Contents**

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



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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