

AhR Bioassay for Environmental Sample Monitoring

The principal application of this Aryl Hydrocarbon Receptor bioassay is to screen processed environmental samples, with the aim of quantifying the aggregate bioactivities, if any, of contaminating biohazardous chemical entities. Contaminants of particular concern include the polycyclic aromatic hydrocarbons (PAHs) and halogenated aromatic hydrocarbons (HAHs) such as dioxin, dioxin-like chemicals, and PCBs. This assay utilizes proprietary mammalian reporter cells expressing the Aryl Hydrocarbon Receptor (AhR; AHR) and a luciferase reporter gene functionally linked to an AhR-responsive promoter. Thus, quantifying changes in luciferase expression in the treated reporter cells provides a sensitive surrogate measure of the changes in AhR activity.

Ordering Information

[Order Online »](#)

NGO-EM06001-0096	96 wells
------------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping

Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name

Aryl Hydrocarbon Receptor

Contents

This Aryl Hydrocarbon Receptor Bioassay for Environmental Sample Monitoring kit is an all-inclusive luciferase reporter assay system that includes, in addition to AhR Reporter Cells, this kit includes are two optimized media for use during cell culture and in diluting the user's test samples, Luciferase Detection Reagent, 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)