

# Zebrafish AR Reporter Assay System

INDIGO's Zebrafish Androgen Receptor Reporter Assay System utilizes proprietary mammalian cells engineered to provide constitutive, high-level expression of Zebrafish (*Danio rerio*) Androgen Receptor (nr3c4), a ligand-dependent transcription factor.

## Ordering Information

[Order Online »](#)

NGO-Z03001-0096	96 wells
NGO-Z03001-0032	3x32 assays

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Shipping                      Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name**                      Androgen Receptor

**Contents**                      This kit product is an all-inclusive assay system that includes, in addition to zAR Reporter Cells, two optimized media for use during cell culture and (optionally) in diluting the test samples, a reference agonist, 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)