

# Validation data for L-Kynurenine

<https://www.invivogen.com/kynurenine>

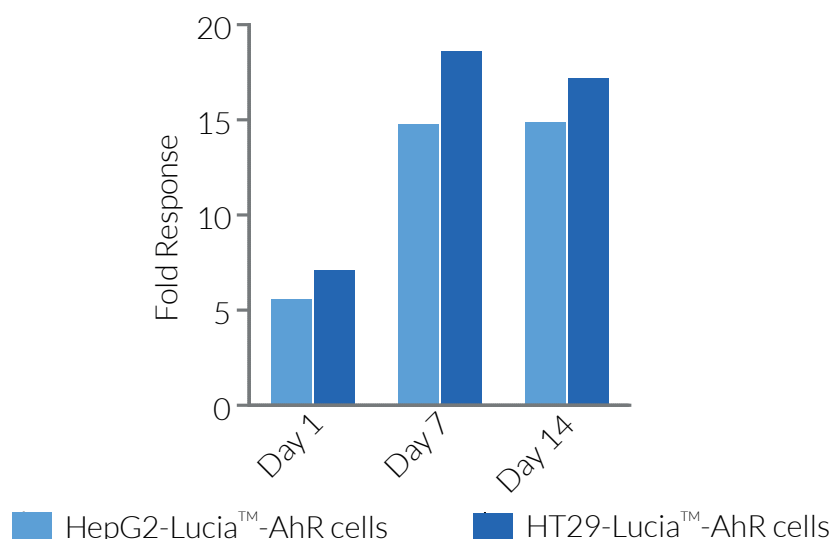
For research use only

Version 19C08-NJ

L-Kynurenine ( $\beta$ -Anthraniloyl-L-Alanine) is an endogenous agonist for the cytosolic aryl hydrocarbon receptor (AhR) that results from tryptophan metabolism. Yet, L-Kynurenine's mode of action on AhR is still unclear. Indeed, L-Kynurenine could be a low-affinity AhR pro-ligand that is slowly converted in high-affinity compounds acting as AhR agonists at subnanomolar concentrations. This conversion occurs independently of enzymes, when fresh crystalline L-Kynurenine is resuspended as a solution, and gives rise to trace-amounts of derivatives named TEACOPs (trace-extended aromatic condensation products).

This increase in L-Kynurenine's potency is detected at least 24h upon resuspension when tested on InvivoGen's HepG2-Lucia<sup>TM</sup>-AhR and HT29-Lucia<sup>TM</sup>-AhR cells. L-Kynurenine's potency increases within days and reaches a plateau between day 7 and day 14 post-resuspension.

## Evaluation of cellular AhR activation by L-Kynurenine



L-Kynurenine induces a time-dependent response in HepG2-Lucia<sup>TM</sup> AhR and HT29-Lucia<sup>TM</sup> AhR cells. The cells were incubated with 33  $\mu$ g/ml (160  $\mu$ M) of L-Kynurenine that had been resuspended and incubated at room temperature for 1, 7 or 14 days. After overnight incubation, the AhR activation was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc<sup>TM</sup>. Data are expressed as fold responses, as compared to non-induced cells. DMSO final concentration impacts cells viability and AhR activation. We recommend to include a control dilution of DMSO in your assay. Using our protocol, a 33  $\mu$ g/ml (160  $\mu$ M) L-Kynurenine concentration includes 1.7% DMSO.

### TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)