

Pollutant Testing & Safety Screening

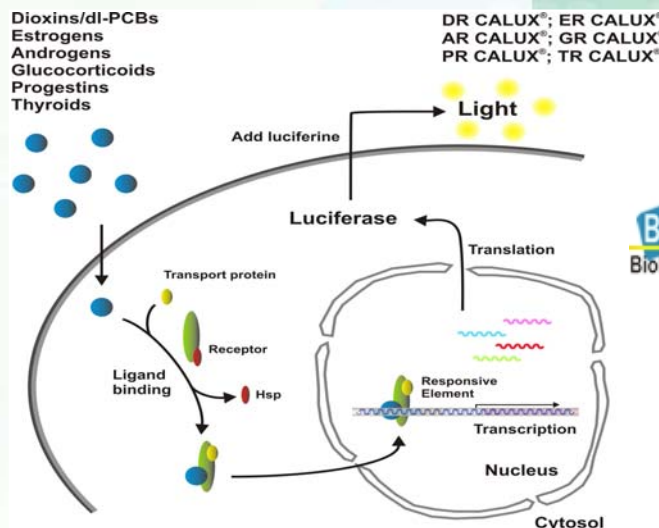
With CALUX[®] Assays and CentroXS³ Microplate Luminometer



The CALUX[®] (**C**hemical **A**ctivated **L**uciferase gene **eX**pression) Assays from BioDetection Systems can be used to detect certain chemicals like dioxins, dioxin-like PCB's or (anti)estrogens and (anti)androgen compounds.

[Learn more about CALUX](#)

The innovative CALUX[®] cells have been tailored so that they produce light in a dose responsive way when exposed to the chemicals. The luciferase activity is quantitated using the BERTHOLD CentroXS3 microplate luminometer



Recommended CentroXS3 model for use with CALUX[®] Assays:

- "Reporter Gene" model (46970-50)
- 2 automatic reagent injectors with JET injection technology (99% precision and 98% accuracy)
- automatic waste pump
- crosstalk: 10⁻⁶
- highest sensitivity (<1 zmol firefly luciferase)



[Learn more about CentroXS3](#)

[Get in touch with Berthold Technologies](#)

www.Berthold.com/Bio