

Extra Sensitivity Centro XS³ LB 960 Microplate Luminometer



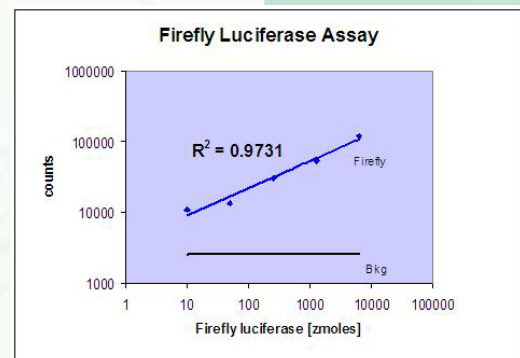
Are you looking for **utmost sensitivity** for your luminescent microplate assays?

You can get detection limits of

- 1 attomol [ATP](#)
- 1 zeptomol [Firefly](#) (~600 molecules)

Have a look at the extra sensitive microplate luminometer [Centro XS³](#).

A new optical design has been implemented to enhance the Centro's detection efficiency even further. In addition this unique system basically eliminates crosstalk to a non-detectable level of 10^{-6} . In addition a most stringent multiplier selection ensures best detection performance.



With the [Centro](#) and the [Centro XS³](#) you can choose ideally suited instruments for

Phagocytosis enzyme kinetics
 Calcium monitoring *reporter gene assays*
ATP determination kinases **caspases**
 toxicity **immuno assays** and more

Consult the [Centro Selection Guide](#) for details.

[Learn more about Centro XS³](#)

[Learn more about Luminescence Applications](#)

[Get in touch with Berthold Technologies](#)

www.Berthold.com/Bio