

SpectraMax Microplate Readers

Comparing luminescence sensitivity of SpectraMax microplate readers using a homogeneous ATP assay

Assay used: ATPlite 1step Luminescence Assay System (PerkinElmer cat. #6016736)

ATP standards ranging from 1 μ M to 1 pM were prepared as a 10-fold serial dilution series in phosphate-buffered saline (PBS). 50 µL of standard was pipetted into the wells of a white 96-well half-area microplate (Corning cat. #3693; n = 4 for each standard). Blank wells containing 50 μ L culture medium only (n = 8) were also prepared. 50 μ L of ATPlite 1step reagent was added to each standard and blank well. Contents of wells were mixed for 2 minutes using a plate shaker. Luminescence was then measured using the microplate readers listed. An integration time of 1 second was used for each reader. Standard curves were plotted using SoftMax® Pro Software (Figure 1).

To assess the sensitivity of each microplate reader, a plate was set up with 32 wells containing PBS only (blank) and four wells containing 50 nM ATP (sample). The lower limit of detection (LLD) for ATP was estimated for each microplate reader using the following formula: LLD = 3*SD_{blank}*Concentration/(RLU_{sample}-RLU_{blank}).

Instruments compared

SpectraMax i3x

SpectraMax® iD

- Onboard luminescence
- Detection cartridge (LUM 96)
- SpectraMax iD5
- SpectraMax L
- SpectraMax M5e
- SpectraMax Mini

ATP LLD

(pM)

59.2

121.3

81.3

36.5

118.1

17.8

Table 1. Lower limits of detection for ATP determined using the ATPlite

Results are shown in Table 1.



Figure 1. ATP standard curves read on SpectraMax microplate readers.

Contact Us

Phone: +1.800.635.5577 Web: www.moleculardevices.com

Email: info@moldev.com Check our website for a current listing of worldwide distributors.

Regional Offices USA and Canada +1.800.635.5577 United Kingdom +44.118.944.8000 00800.665.32860 Europe* China +86.4008203586

Taiwan/Hong Kong +886.2.2656.7585 Japan +81.3.6362.9109 +82.2.3471.9531 South Korea +91.73.8661.1198

1step assay with SpectraMax microplate readers.

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom

Instrument

SpectraMax i3x,

SpectraMax iD5

SpectraMax M5e

SpectraMax Mini

India

SpectraMax L

SpectraMax i3x, luminescence

detection cartridge LUM96

onboard luminescence

The trademarks used herein are the property of Molecular Devices, LLC or their respective owners. Specifications subject to change without notice. Patents: www.moleculardevices.com/productpatents FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

ATP LLD

(moles)

5.9 x 10⁻¹⁵

 1.2×10^{-14}

8.1 x 10⁻¹⁵

 3.6×10^{-15}

1.2 × 10⁻¹⁴

1.8 x 10⁻¹⁵