

# Can it even read Chinese?



## After reading in different modes, she asks herself if there is anything the POLARstar Omega can't read

The new POLARstar Omega can't read Chinese, but it is the first multidetection microplate reader to offer high speed, full spectrum absorbance measurements. For the greatest flexibility in microplate-based assays, four additional detection modes are also available: fluorescence intensity, time-resolved fluorescence, fluorescence polarization and high performance luminescence. The new spectrometer technology allows full UV/Vis spectrum analysis from

220 to 850 nm at a resolution of 1 nm. Read times faster than 1 second per well are possible. Simultaneous dual emission detection in fluorescence and luminescence modes complete the POLARstar Omega reader. Because of our continuous innovations, BMG LABTECH has become a world leader in microplate reading systems, offering a broad portfolio of solutions for your life science research. With the new POLARstar Omega you are never lost in translation.

Microplate readers for all your assay needs



FLUOstar

PHERAstar

NOVOstar

NEPHELOstar

Stacker

**BMG LABTECH**

[www.bmglabtech.com](http://www.bmglabtech.com)