

## SPECTROstar® Nano - Technical specifications

Due to the modularity of BMG LABTECH's instruments, all or combinations of the features below can be installed at purchase or upgraded at any time. Please contact your local representative for more details or a quote.

<b>Detection modes</b>	UV/vis absorbance spectrum
<b>Measurement modes</b>	Endpoint and kinetic measurements Well scanning
<b>Microplate formats</b>	6- to 1536-well plates, user-definable
<b>Microplate carrier</b>	Robot compatible
<b>Cuvette port</b>	Cuvette port for cuvettes with 10 mm path length Micro cuvettes, Traycell compatible Beam height 8.5 mm
<b>Light source</b>	High energy xenon flash lamp
<b>Detector</b>	Spectrometer with CCD
<b>Spectral range</b>	220 - 1000 nm
<b>Sensitivity</b>	Selectable spectral resolution: 1, 2, 5, 10 nm OD range: 0 to 4 OD Accuracy: < 1% at 2 OD Precision: < 0.5% at 1 OD and < 0.8% at 2 OD
<b>Read times</b>	Full spectrum from 220 to 1000 nm in less than 1 sec/well
<b>Shaking</b>	Linear, orbital, and double orbital User-definable time and speed
<b>Incubation</b>	+3°C above ambient up to 45°C for microplate and cuvette
<b>Gas vent</b>	System to inject an atmosphere or to pull a vacuum into the reader
<b>Software</b>	Multi-user software package including Reader Control and MARS data analysis software, FDA 21 CFR part 11 compliant
<b>Computer interface</b>	USB 2.0 compatible to USB 1.1
<b>Dimensions</b>	Width: 36 cm, depth: 50 cm, height: 16 cm Weight: 10 kg
<b>Accessories</b>	
<b>LVis Plate</b>	Microplate designed to measure 16 low-volume (2 µL) samples and standard cuvettes. Incorporates NIST-traceable filters and holmium oxide standards for instrument performance testing. Sensitivity: 2 ng/µL dsDNA
<b>Barcode reader</b>	External manual barcode reader
<b>THERMOstar</b>	Microplate incubator and shaker

Specifications are subject to change.  
Limit of detection (sensitivity) was calculated according to the IUPAC standard:  $3 \times \text{SD}_{\text{blank}} / \text{slope}$   
SPECTROstar is a registered trademark of BMG LABTECH GmbH.  
© 2016 All rights reserved. All logos and trademarks are the property of BMG LABTECH.