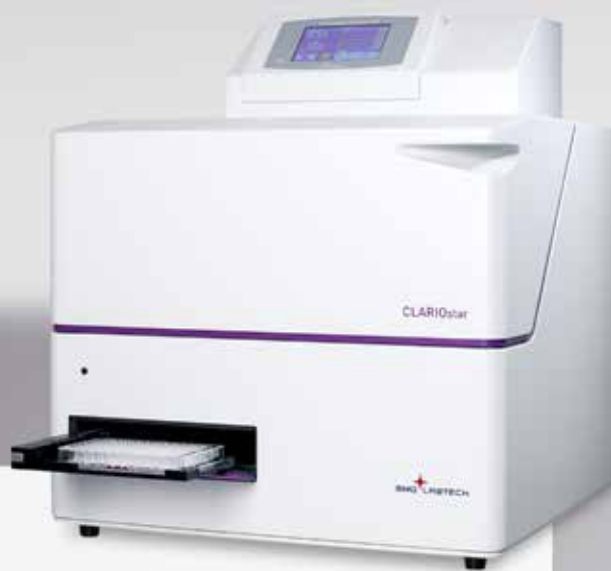


# CLARIOstar®

with Atmospheric  
Control Unit (ACU)



## Independent gas control for happy cells

The Atmospheric Control Unit (ACU) module allows independent, simultaneous regulation of O<sub>2</sub> and CO<sub>2</sub> within the microplate reader chamber. From standard cell growth to hypoxic or cytotoxicity assays, the ACU provides the optimal environment for any live cell-based assay.

### Features:

- O<sub>2</sub> and CO<sub>2</sub> regulation
- Gas value trackability in MARS data analysis software
- O<sub>2</sub> and CO<sub>2</sub> gas ramps
- Altitude correction for accurate gas regulation
- Intuitive interface with touchscreen, up to 10 gas presets and gas concentration curve display



## Technical Specifications

$O_2$ and $CO_2$ control	Range: 0.1 – 20 % Control: $\pm 0.1$ % Sensor: Low drift, long lifetime
--------------------------	---

All active components are integrated inside the ACU for highest reliability, independence of the microplate reader and easy maintenance. In combination with temperature control, three different shaking options, bottom reading detection, and Z-height focus adjustment, the ACU provides an ideal 'walk-away' solution for any long-term cell-based assay.

## Applications

- Proliferation assays and cell viability
- Migration and invasion studies
- Hypoxia and Ischemia/reperfusion
- Angiogenesis
- Cytotoxicity studies
- Viral uptake