

Atmospheric Control Unit

for Omega series and NEPHELOstar^{Plus}



Independent gas control for happy cells

The Atmospheric Control Unit (ACU) allows independent, simultaneous regulation of O₂ and CO₂ within the microplate reader.

From standard cell growth to hypoxic or cytotoxicity assays, the ACU provides the optimal environment for any live cell-based assay.

Features:

- Independent O₂ and CO₂ regulation
- Up to 10 user-definable gas presets
- Touchscreen with intuitive interface
- Gas concentration curve display
- Altitude correction for accurate gas regulation
- Compatible with SPECTROstar[®], LUMIstar[®], FLUOstar[®] and POLARstar[®] Omega, and NEPHELOstar^{® Plus}

Technical Specifications

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

O₂ and CO₂ control	Range: 0.1 – 20 % Control: ± 0.1 % Sensor: Low drift, long lifetime
---	---

Applications

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

