

# Luminex® XYP™



## Specifications

### Heater Operating Range

35°C to 55°C (95° to 131° F)  
with tolerance 0°C to +2°C.

### Physical Dimensions

44 cm (W) x 61 cm (D) x 8 cm (H) (approx.)  
17.25" (W) x 24" (D) x 3" (H) (approx.)  
Weight: 14.5 kg (32 lbs.)

### Electrical

Power: 100-240 VAC, 2.25 Amps, 47-63 Hz

### Communications Interface

RS--232

### Operating Environment

Operating temperature 15°C to 30°C (59° to 86° F)  
10% - 90% relative humidity non condensing  
Altitude: up to 2400 m (7874ft) above mean sea level or below.

### 96 Well Microtiter Plates

Compatible with standard plates no thicker than 0.75 inches (19mm)

### Order Information

Luminex XYP  
Includes accessories  
Product No. CN-X004-01  
To order, call: 866-401-5450

For technical support, call: 877-785-2323 (U.S. and Canada) or +1 512-381-4397 (International)

Fax: 512-219-5114

E-mail: [support@luminexcorp.com](mailto:support@luminexcorp.com)

**Luminex Corporation**  
12212 Technology Blvd.  
Austin, Texas 78727  
[www.luminexcorp.com](http://www.luminexcorp.com)