

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 2000 m (6600ft) 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-X001-001

To order, call: 888-219-8020

For technical support, call: 877-785-2323 (U.S. and Canada) or +1 512-381-4397 (International)
Fax: 512-219-5195 E-mail: support@luminexcorp.com

Luminex Corporation
12212 Technology Blvd.
Austin, Texas 78727
www.luminexcorp.com