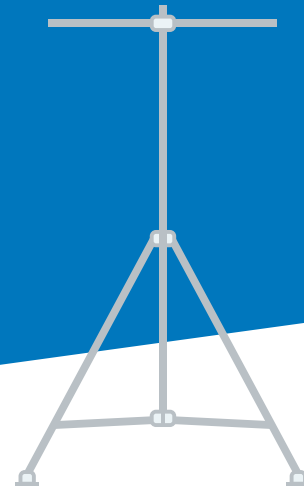


Tripod and Mounting Hardware

A Solid Foundation for Your Research



Protect your investment with the rugged heavy duty adjustable tripod. This sturdy instrument platform provides strong support for eddy covariance systems, meteorological sensors, and enclosures in long-term field deployments.

Designed to withstand extreme outdoor conditions, including high winds, the tripod extends to a height of up to 4 meters and can withstand hurricane-force winds of up to 100 mph with a full load of mounted instruments.

To learn more about the tripod and mounting options for your eddy covariance system visit [licor.com/comms_power](https://www.licor.com/comms_power).

Key Features

- Confirmed by certified, independent engineers to withstand 100 mph winds with a full load of instruments
- Fully adjustable up to 4 meters
- Includes all necessary guy wires and a lightning rod
- Includes carrying case to aid with portability

Tripod and Mounting Hardware Specifications

Tripod

Load Capacity: 135 kg (297 lbs)

Maximum Load Area: 2.19 m² (23.58 ft²)

Base Diameter: 2.54 m (8 ft 4 in)

Mast Outer Diameter: 6.03 cm (2.375 in)

Weight: 50 kg (110 lbs)

Maximum Slope: 13.0°

Maximum Wind Gust (3 seconds): 161 kph (100 mph)*

* When anchored in recommended soil conditions.

Upper Mast

Materials: 6061-T6 Aluminum Pipe (2-inch Iron Pipe Size, Schedule 40)

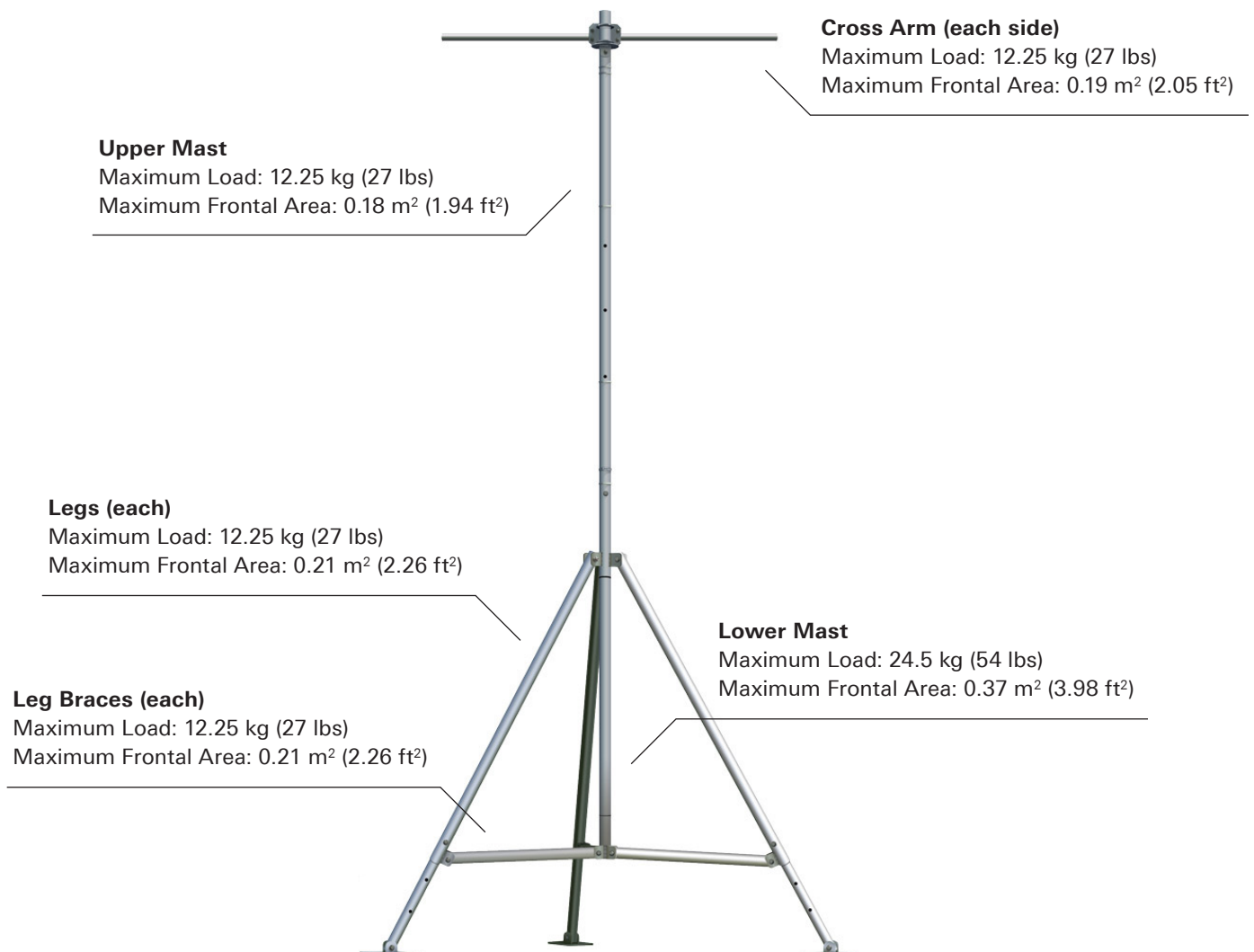
Outer Diameter: 6 cm (2.375 inches)

Cross Arm

Materials: 6061-T6 Aluminum Pipe (1-inch Iron Pipe Size, Schedule 80)

Outer Diameter: 3.3 cm (1.315 inches)

Length: 1.5 m (59 inches)



Specifications subject to change without notice.