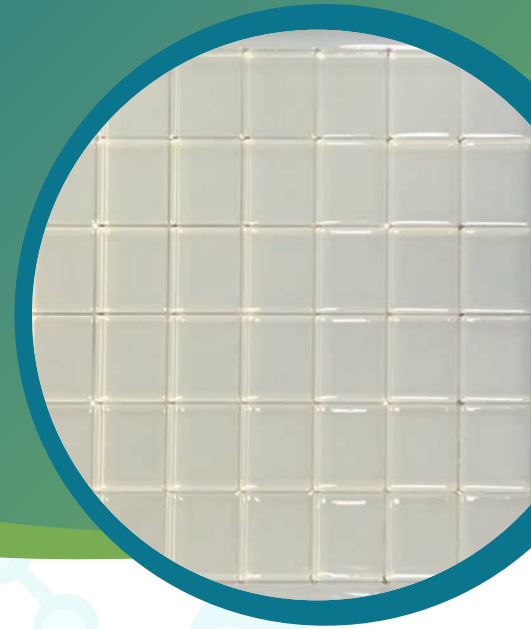


**DATA SHEET**

# Vented Q-Tray with Cover and 48 Well Divider

Catalog Part Number: X6029

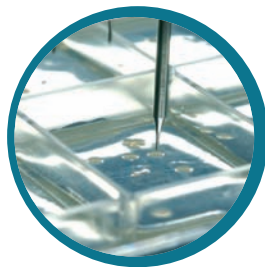
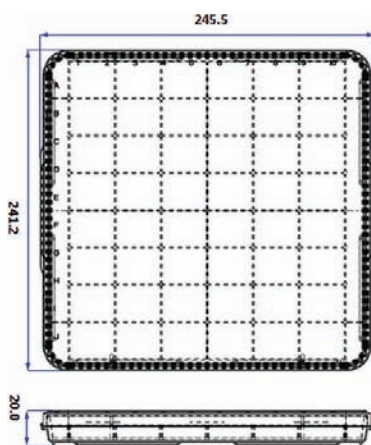


## Material(s)

Optically Clear Polystyrene for Q-tray Dish/Cover/Divider

## Assembled product dimensions

- 245.1 mm (L) x 241mm (W) x 20mm (H)
- Tolerances:  $\pm 0.4$ mm



## Sterilization

The product has been Irradiated and dosimetrically released in accordance with ISO 11137 – *Sterilization of Health Care Products*

## Performance testing

Each manufactured lot is sampled and tested in accordance with Standard Operation Procedures.

## Visual attributes

- Visual and tool-aided inspection of the product appearance, dimensions, and optical clarities
- Inspection for compatibility of the product with QPix automated colony picker instruments

## Packaging

- Inspections for correct labeling, configuration, and package seal integrity

## Hot agar tolerance

- Up to 68 °C / 154.4 °F

## Contact Us

Phone: +1.800.635.5577  
 Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
 Email: [info@moldev.com](mailto:info@moldev.com)  
 Check our website for a current listing of worldwide distributors.

## Regional Offices

USA and Canada	+1.800.635.5577	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.3.6362.9109
Europe*	00800.665.32860	South Korea	+82.2.3471.9531
China	+86.4008203586	India	+91.73.8661.1198

\*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom