

QReps

Disposable Polypropylene Replicators

Molecular Devices QReps are a range of polypropylene handheld replicators that can be used for a range of applications including the manual replication of bacterial clone libraries and the manual printing of microarrays for screening via hybridization.

The disposable plastic QReps are made from polypropylene and are supplied sterile, eliminating the need for flaming or ethanol washing prior to use. A choice of 96- or 384-pin versions are available in different pin lengths and thickness for maximum user choice and application flexibility. The 10 mm pin length QReps are compatible with Molecular Devices low profile plates while the 15 mm pin length is best suited to standard deep-well plates.

All QReps and hand held replicators can be used with the Molecular Devices microplates and deep-well plates.

Technical information

Material: polypropylene.

Sterility: supplied sterile (gamma irradiated)

Chemical resistance: High (see polypropylene chemical resistance chart)

Polypropylene chemical resistance:

Resistant to the following chemicals at room temperature.

- Acetic acid
- Acetone
- Acetonitrile
- Amino Acids
- Ammonia
- Ammonium Acetate
- Aniline (50% wt)
- Benzene
- Butanol
- Calcium Hydroxide
- Carbazole
- Chloroacetophenone
- Chloroform
- Cyclohexanol
- Cyclohexanone
- Diethylene Glycol
- Dimethylformamide (DMF)
- Dimethylsulfoxide (DMSO)
- Dipropylene Glycol
- Ethanol
- Ethyl Alcohol
- Ethyl Acetate
- Ethylene Glycol
- Fluorides
- Formaldehyde (30%)
- Formic Acid
- Hydrochloric Acid (30%)
- Hydrofluoric Acid
- Hydrobromic Acid (35%)
- Isopropyl Alcohol
- Methanol
- Methyl Acetate
- Methylene Chloride
- Phosphoric Acid (50%)
- Potassium Hydroxide (10%)
- Propanol
- Sodium Acetate
- Sodium Chloride
- Sodium Hydroxide (15%)
- Sodium Nitrate (10%)
- Sulphuric Acid
- Tetrahydrofuran (THF)
- Toluene
- Triethelene Glycol

Source: <http://www.vp-scientific.com/polypropylene%20properties.htm>

Ordering information

Catalog No.	Description	Pack size
X5050	QRep 384 Pin Replicators: 10 mm length, 1 mm diameter	128
X5051	QRep 96 Pin Replicators: 10 mm length, 1 mm diameter	128
X5052	QRep 96 Pin Replicators: 10 mm length, 2 mm diameter	128
X5053	QRep 384 Pin Replicators: 15 mm length, 1 mm diameter	112
X5054	QRep 96 Pin Replicators: 15 mm length, 1 mm diameter	112

Contact Us

Phone: +1-800-635-5577
 Web: www.moleculardevices.com
 Email: info@moldev.com
 Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada +1-800-635-5577
 Brazil +55-11-3616-6607
 China (Beijing) +86-10-6410-8669
 China (Shanghai) +86-21-3372-1088
 Germany 00800-665-32860

Japan (Osaka) +81-6-7174-8831
 Japan (Tokyo) +81-3-6362-5260
 South Korea +82-2-3471-9531
 United Kingdom +44-118-944-8000