

## Thermo Scientific HERAfreeze HDE Series Ultra-Low Freezers, 600-Box Capacity, 230V, 50Hz

Asheville, North Carolina, United States

SKU HDE60086LV



### Manufacturing

Manufacturing Impact Reduction	3.0
Renewable Energy Use	No
Responsible Chemical Management	1.0
Shipping Impact	9.3
Product Content	5.0
Packaging Content	5.0

### User Impact

Energy Consumption (kWh/day)	12.7
Water Consumption (liters/day)	N/A
Product Lifetime	5.0

### End of Life

Packaging	5.3
Product	1.0

### Innovation

Innovative Practices	-1.0
----------------------	------

**Environmental Impact Factor: 46.3**

Label Valid Through:

June 2023

