

Thermo Scientific TDE Series Ultra-Low Freezers, 600-Box Capacity, 230V, 50Hz

Asheville, North Carolina, United States

SKU TDE60086FV

**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

