

## Thermo Scientific TDE Series Ultra-Low Freezers, 600-Box Capacity, 115V, 60Hz

Asheville, North Carolina, United States **SKU** TDE60086FA

222	Environmental Impact Scale	
<b>←</b>	Decreasing Environmental Impa	ct 10
Manuf	facturing —	
Manufacturing Impact Reduction		3.0
Renewable Energy Use		No
Responsible Chemical Management		1.0
Shipping Impact		1.0
Product Content		5.0
Packaging Content		5.0
User I	mpact —	
Energy Consumption (kWh/day)		12.7
Water Consumption (gallons/day)		N/A
Produc	t Lifetime	5.0
End o	f Life —	
Packaging		5.3
Produc	t	1.0
Innova	ation —	
Innovative Practices		-1.0
Enviro	onmental Impact Factor:	38.0
Label Va	alid Through:	June 2023



act.mygreenlab.org