

Introducing the ACT Ecolabel by My Green Lab

What Is the ACT Ecolabel?

The ACT Ecolabel program by My Green Lab aims to address the need of both scientists and procurement specialists for clear, third-party verified information about the environmental impact of laboratory products.



The ACT Ecolabel, which stands for Accountability, Consistency, and Transparency (ACT), is a standardized measure of a product's environmental impact, assisting laboratories in making informed, sustainably conscious choices when comparing products from different manufacturers.

Key Aspects Addressed by the ACT Ecolabel:



Manufacturing: The product's construction, content, and the manufacturer's environmental policies.



Packaging: The materials used, and the waste generated by the packaging process.



Energy and Water Use: The product's efficiency throughout its operational lifespan (where applicable).



End-of-life: The product's disposal, recyclability, and reusability options.



Shipping: Impact of shipping the product from the site of manufacturing to end customer.



Innovation: Advancing sustainability practices to enhance the product's environmental performance and minimize its ecological footprint.

How Does it Work?

Products with the ACT Ecolabel have undergone independent auditing by SMS Collaborative, LLC (SMSC) to ensure that the provided information is accurate and impartial. My Green Lab makes the ACT Ecolabels available in a publicly accessible database:

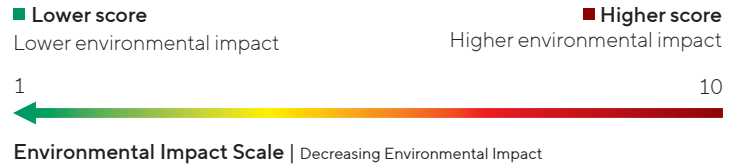
Act - ACT Database (mygreenlab.org)

Why ACTing Matters?

At Sartorius, we share your commitment to sustainability and are dedicated to maintaining transparency in our manufacturing processes and product content. Our collaboration with the My Green Lab program and the adoption of the ACT Ecolabel are concrete steps we have taken to ensure open and clear communication in laboratory procurement. The time to ACT is now!

How to Read the ACT Ecolabel

Each product category is evaluated on a scale from 1 to 10, with a lower score indicating a smaller environmental footprint. The criteria for the ACT Ecolabel are formally referred to as the Environmental Impact Factor (EIF). Products that achieve a lower EIF score have demonstrated a reduction in their environmental impact through various means, such as using renewable energy sources, minimizing waste, utilizing recyclable materials, and implementing efficient manufacturing processes.



This category is either Yes or No.



ACT. The Environmental Impact Factor Label		US
Product Name		
Product Location SKU 0000		
1 ← 10 		
Environmental Impact Scale		
Decreasing Environmental Impact		
Manufacturing		
Manufacturing Impact Reduction	3	
Renewable Energy Use	Yes	
Responsible Chemical Management	5	
Shipping Impact	9	
Product Content	1	
Packaging Content	5	
User Impact		
Energy Consumption (kWh/day)	2.5	
Water Consumption (gallons/day)	13.1	
Product Lifetime	4	
End of Life		
Packaging	5	
Product	1	
Innovation		
Innovative Practises	-1	
Environmental Impact Factor	47,6	
Label Valid Through:	January 2021	
act.mygreenlab.org		



Regional labels reflect the differences in Shipping and End-of-Life impacts for each region (US, EU, UK) where the product is sold. They also highlight how energy usage may vary across markets and report water usage in units specific to each market (gallons per day or liters per day) for equipment.



These values are graded on a scale of 1-10, with 1 indicating the lowest environmental impact and 10 indicating the highest environmental impact.



These values represent actual daily consumption and apply only to equipment/instruments.



The sum of all values equals the Environmental Impact Factor. A lower number indicates a lower overall environmental impact.

ACT Ecolabels are valid for two years from the date of issue.



Summary

The ACT Ecolabel serves as a tool for those looking to make more sustainable choices in their laboratories. It simplifies the decision-making process by providing clear, standardized information on the environmental impact of laboratory products. Choosing products with lower environmental impact scores allows labs to contribute to a collective effort to minimize the ecological footprint of scientific research and laboratory work. This transparency enables easy comparison of products and informed decision-making that aligns with personal and laboratory sustainability goals, ultimately facilitating a more straightforward and responsible procurement process.