

# World Hunger Facts

visit [bio-rad.com/worldhunger](http://bio-rad.com/worldhunger)

Over 800 million people worldwide are chronically undernourished



One in 10 experiences chronic hunger.



60% of the world's hungry are women



Over 40 million Americans live in food-insecure homes.



**United Nations Sustainable Development Goal:**  
End hunger and achieve food security with better nutrition.



**How?**  
Sustainable systems of production, distribution, and consumption of food.

## Environmental

Develop adaptations to climate change, extreme weather, drought, and flooding. Maintain genetic diversity of seeds, cultivated plants, and farmed/domesticated/wild animal species. Improve soil quality.



## Social

Ensure access by all people to safe, nutritious, and sufficient food year-round.



## Economic

Implement agricultural systems for increased productivity and production for:

- Small-scale food producers
- Indigenous people
- Family farmers

Limit trade restrictions in global agricultural markets.

Bring engineering into your classroom as students tackle the global issue of world hunger. Visit **bio-rad.com/worldhunger** for more information.

### Ordering Information

<i>Catalog #</i>	<i>Description</i>
17005278EDU	<b>Engineering Solutions for Global Health Kit</b>



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00  
**Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600  
**Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30  
**Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580  
**Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300  
**United Kingdom** 44 01923 47 1301

