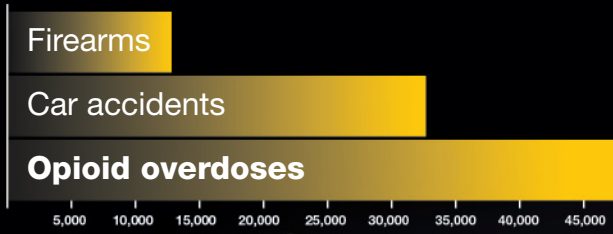


The Opioid Epidemic in the U.S.

bio-rad.com/opioidinfo

Opioid drug overdoses kill more people than car accidents and firearms combined.



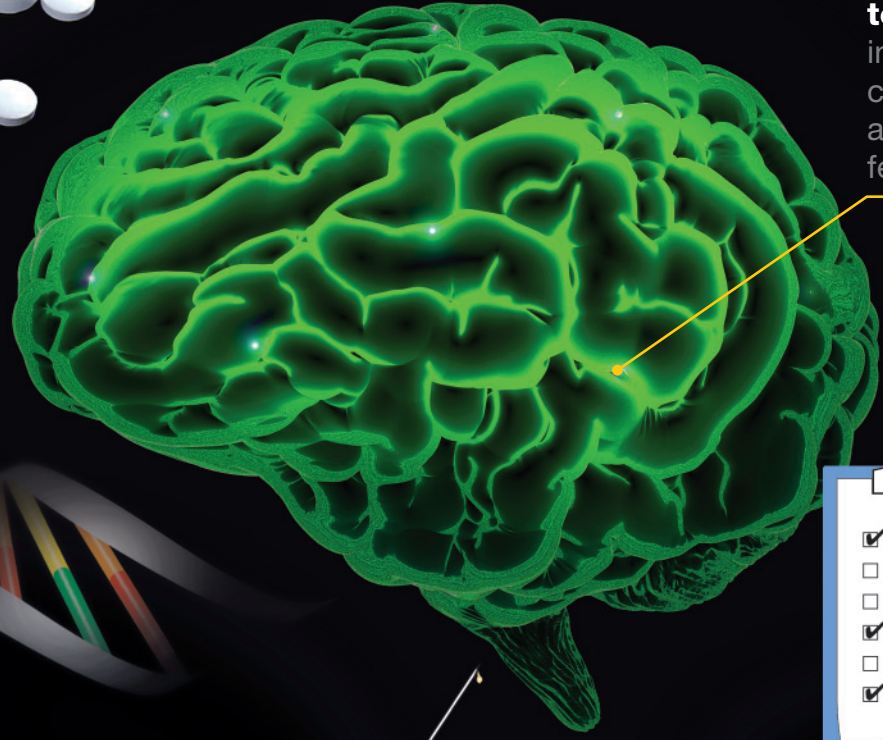
The opioid crisis

costs the U.S. **\$75 BILLION** annually.



17%

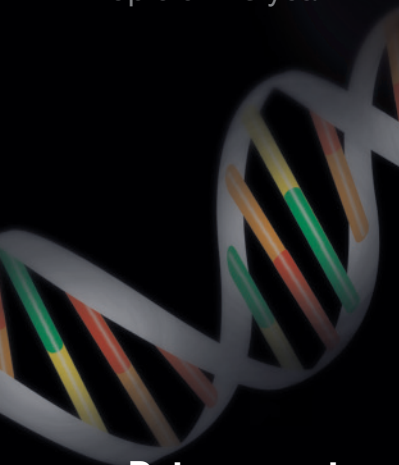
of Americans had at least one prescription opioid this year.



Opioids bind to receptors in the brain, spinal cord, and other areas and reduce feelings of pain.

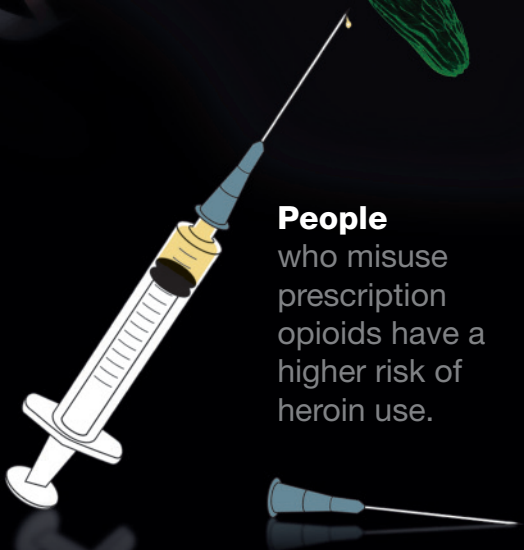
-
- Medication
 - Hospital Stay
 - Residential Stay
 - Therapy
 - MAT
 - Counseling

Treatment for opioid abuse includes medication, residential/hospital stays, medication-assisted therapy (MAT), and counseling.



Data suggest complex genetic and environmental influences on opioid dependence.

People who misuse prescription opioids have a higher risk of heroin use.



Bring data and research sciences into your classroom as students tackle the global issue of opioid dependence. Visit bio-rad.com/opioid for more information.

Ordering Information

<i>Catalog #</i>	<i>Description</i>
17005316EDU	Science of Opioid Dependence Kit
17005313EDU	Science of Opioid Dependence Kit plus UView Electrophoresis Reagent Pack
17005297EDU	Science of Opioid Dependence Kit plus Fast Blast Electrophoresis Reagent Pack

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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

