

Biofuel Enzyme Reactions Kit

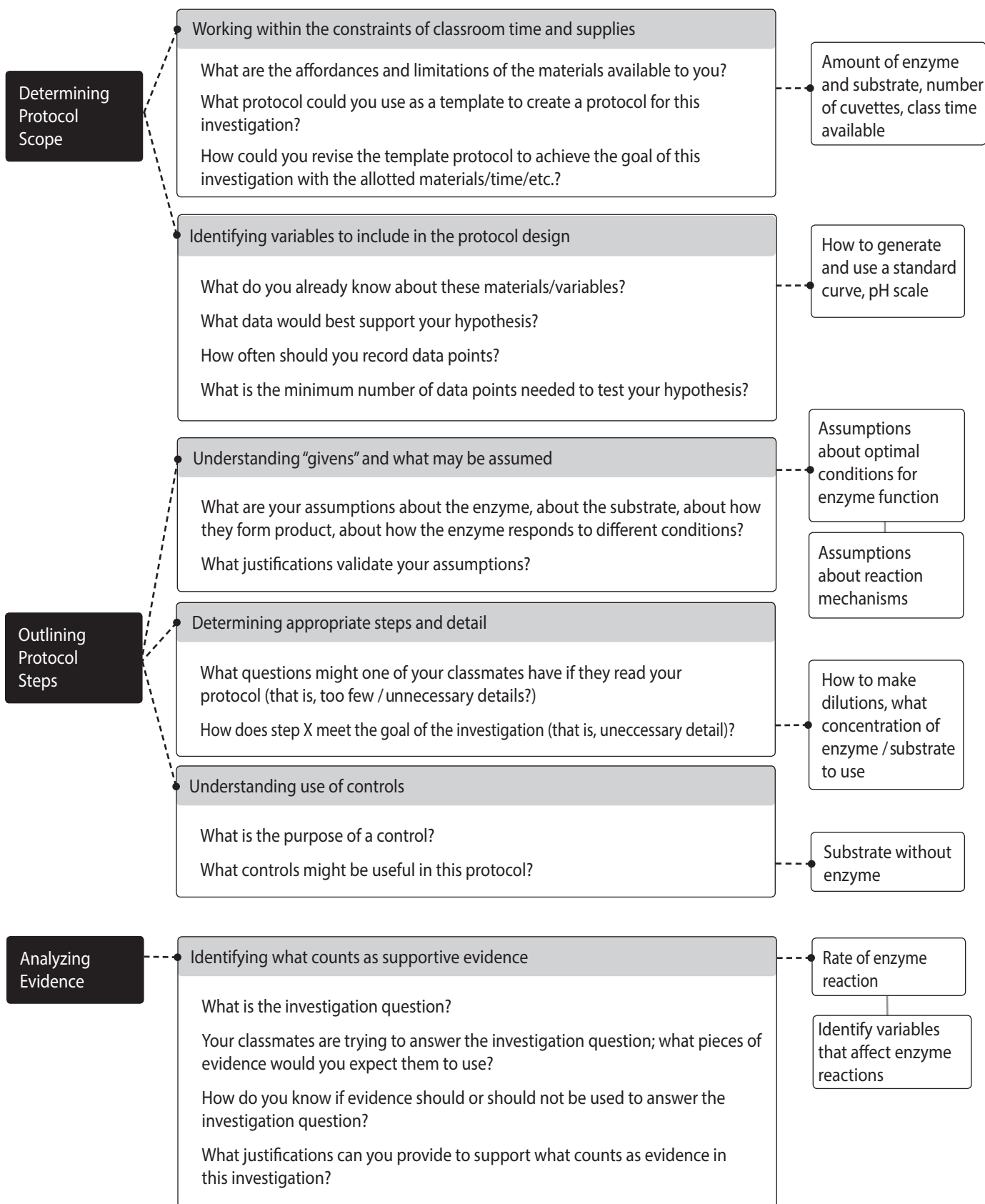
A ThINQ![™] Investigation

Teacher Model Process

Guide to Student Inquiry and Investigation

This table is designed to highlight specific steps during protocol design where students may require additional support. As students design their protocols you may find it useful to support their thinking and writing by using the questions and prompts below. This table can be used in conjunction with the Experimental Planning and Design Worksheet (bio-rad.com/doc/biofuelAPresources) as a formative or summative assessment tool and during class time to support students in the protocol design process.

Inquiry Lesson Step	Suggested Questions and Prompts to Support Protocol Design	Kit Specific Applications
Making Observations	<p>Making observations that lead to an investigation question</p> <p>What did you notice during Investigation 1 when using mushroom extract to determine reaction rate?</p> <p>How might different conditions affect the rate of the enzyme reaction?</p> <p>What kinds of observations would you need to make in order to answer your questions?</p> <p>Describe the phenomenon that you observed in Investigation 1.</p>	<p>Determining reaction rate using enzyme extracted from mushroom, comparing enzyme activity from different mushrooms</p>
Defining the Purpose of the Investigation	<p>Clarifying the purpose of the investigation</p> <p>What was the purpose of Investigation 1?</p> <p>How does the purpose of Investigation 1 differ from the purpose of this investigation?</p> <p>What small changes could you make to the protocol in Investigation 1 to meet the purpose of this investigation?</p> <p>What steps from the protocol in Investigation 1 can you use to design this investigation?</p>	<p>Comparing different conditions of temperature, pH, enzyme concentration, and substrate concentration</p>
Hypothesis Formation	<p>Clarifying goals for the investigation</p> <p>Can you explain in your own words what the investigation question is asking?</p> <p>What do you already know about how enzymes work?</p> <p>How does your knowledge about enzymes help you describe the phenomena that you observed in Investigation 1?</p> <p>Knowing this, how would you modify your model to describe the phenomena that you observe in this investigation?</p> <p>What evidence would you need in order to answer the investigation question?</p>	<p>Understanding how enzymes function under different conditions (e.g., some denature at high temperatures)</p>





**Bio-Rad
Laboratories, Inc.**

*Life Science
Group*

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00
France 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 972 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 63617000 **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 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026674 55 05
Taiwan 886 2 2578 7189 **Thailand** 66 662 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

