



Unit 1: The Relationship between Genes and Proteins

THE RELATIONSHIP BETWEEN GENES AND PROTEINS: TRANSCRIPTION AND TRANSLATION

This unit introduces students to the processes of transcription and translation. The unit includes an accompanying laboratory *Effects of antibiotics on bacterial growth and protein synthesis* to allow students to explore lecture concepts and principles with hands-on exercises.

Lecture Presentation



The Relationship between Genes and Proteins: Transcription and Translation

(1 MB)

Laboratory Protocols



Instructor's Lab Manual

(206 KB)



Student's Lab Manual

(87 KB)

Instructor's Guide



Instructor's Guide for Unit 1

(152 KB)

Animations

Transcription Animation (PC)

(4 MB)

Transcription Animation (Mac)

(25 MB)

Translation Animation (PC)

(6 MB)

Translation Animation (Mac)

(1 MB)

More Transcription and Translation Resources

[All your ribosomes belong to us.](#) This Science Blog does a great job giving perspective to the role of the ribosome in the biology of the cell and making the ribosome exciting.