

# SAMPLE SUBMISSION FORM

Date:

Sender:

Company/Institution:

Contact information:

*(email and or telephone number)*



Date of sample collection	Sample type	Volume	Quantity	Sample ID	Protocol Number / Study Identification / Job ID	Comment / brief protocol description