



BIO-RAD LABORATORIES, INC. POLITICAL ACTIVITY POLICY

I. Introduction

Bio-Rad's mission is to advance science and save lives, together. Bio-Rad is committed to pursuing this mission with integrity, which means conducting ourselves in a professional and ethical manner and complying with all applicable laws and regulations in all that we do. Bio-Rad values diversity and inclusion in our workforce and we encourage our employees to take an active role in community and civic organizations and to make a difference in the issues that matter to them, whatever they might be. Because we value and support a diverse workforce with diverse viewpoints, we believe it is important that Bio-Rad focus on our mission and not, as a company, become involved in political activities. Accordingly, Bio-Rad has a long-standing policy prohibiting the Company, and employees acting in their official capacity, from supporting any political candidate or participating in any political cause or concern.

II. Prohibition on Corporate Political Activity

It is Bio-Rad's policy that the Company's time, funds, or other resources shall not be used, directly or indirectly, to advance or support any political cause, candidate, campaign, or party. Employees should never suggest that they speak for Bio-Rad on any political issue or for or against any political candidate or referendum, and employees shall not make political donations or contributions under Bio-Rad's name. This prohibition includes contributions to political committees (including "527 organizations" such as political action committees and ballot measure committees), to influence the outcome of ballot measures, to trade organizations with the express purpose of direct or indirect political participation or election-related activity, and to tax-exempt groups, such as 501(c)(4) social welfare organizations that may use such funds for political or election-related purposes. Bio-Rad itself does not sponsor a political action committee (PAC) and does not contribute to PACs. While employees are encouraged to get involved and make a difference in the issues that matter to them, when participating in the political process, they must make sure to do so only on their personal time and using only their personal resources.

In accordance with our Global Anti-Bribery & Anti-Corruption Compliance Policy, Company employees may not use expense accounts to pay for any personal political contributions or seek any form of reimbursement from Bio-Rad for such contributions. Furthermore, a Company employee may never make a political contribution, even in his or her own name, to obtain or retain business or to gain an improper business advantage.

Bio-Rad values trade associations and the Company is an active participant in certain select industry groups. Trade associations provide important services to their members, including monitoring legislative and regulatory agendas, providing data and analysis, and educating lawmakers on important scientific advances. Trade associations in which Bio-Rad participates do not use member dues to advance political causes or candidates.