

## Quanterix Board Diversity Matrix as of February 1, 2023

Total Number of Directors	<b>9</b>			
	<b>Female</b>	<b>Male</b>	<b>Non-Binary</b>	<b>Did Not Disclose Gender</b>
<b>Part I: Gender Identity</b>				
Directors	<b>3</b>	<b>5</b>	-	<b>1</b>
<b>Part II: Demographic Background</b>				
African American or Black	-	-	-	-
Alaskan Native or Native American	-	-	-	-
Asian	-	-	-	-
Hispanic or Latinx	-	-	-	-
Native Hawaiian or Pacific Islander	-	-	-	-
White	<b>3</b>	<b>5</b>	-	-
Two or More Races or Ethnicities	-	-	-	-
LGBTQ+	<b>1</b>			
Did Not Disclose Demographic Background	<b>1</b>			