

PUBBLICAZIONE 2020 TRASPARENZA BIO-RAD

I soci devono documentare e rendere pubblici ogni anno attraverso un apposito Modello di Trasparenza che costituisce parte integrante del presente Codice , i trasferimenti di valore effettuati direttamente o indirettamente ai Professionisti del Settore Sanitario, alle Organizzazioni Sanitarie e alle Terze Parti.

	Full Name	HCPs: City of Principal Practice HCOs: City where registered	Country of Principal Practice	Principal Practice Address	Donations	Scholarship	Contribution to costs of Events			Contribution to costs of Training			Costs for the participation in training, educational and promotional events on company's products organised by the Members	Fee for service and consultancy		Transfers of Value Research & Development	Total	
							Sponsorship agreements with HCOs/Third Parties appointed by HCOs to manage an event	Registration Fees	Travel & Accommodation	Sponsorship agreements with HCOs/Third Parties appointed by HCOs to manage an event	Registration Fees	Travel & Accommodation		Fees	Related expenses agreed in the fee for service or consultancy contract, including travel e accommodation relevant to the contract			
HCPs (Healthcare Professionals)	INDIVIDUAL NAMED DISCLOSURE - one line per HCP (i.e. all transfers of value during a year for an individual HCP will be summed up: Itemization should be available for the individual Recipient or public authorities' consultation only, as appropriate)																	
					NA	NA	NA*	NA*	NA*								NA	
					NA	NA	NA*	NA*	NA*								NA	
					NA	NA	NA*	NA*	NA*								NA	
					NA	NA	NA*	NA*	NA*								NA	
	AGGREGATE DISCLOSURE - per HCPs																	
	Aggregate amount attributable to transfers of value to such Recipients				NA	NA	NA*	NA*	NA*								NA	
	Number of Recipients in aggregate disclosure				NA	NA	NA*	NA*	NA*								NA	
	% of the number of Recipients included in the aggregate disclosure in the total number of Recipients disclosed				NA	NA	NA*	NA*	NA*								NA	
	INDIVIDUAL NAMED DISCLOSURE - one line per HCOs (i.e. all transfers of value during a year for an individual HCOs will be summed up: Itemization should be available for the individual Recipient or public authorities' consultation only, as appropriate)																	
Healthcare Third Parties	Selene S.r.l. – Eventi e Congressi	Torino	Italy	Via G. MEDICI 23, 10143 Torino	Only HCO	NA	3,500			NA	NA	NA	NA				NA	
	SIMTIPTRO Srl	Milano	Italy	Via desiderio 21, 20131 Milana	Only HCO	NA	5,500			NA	NA	NA	NA				NA	
	Innovaformez Srl	Catanzaro	Italy	Via Graddoni Ospedale Civile, 11, 88100 Catanzaro	Only HCO	NA	4,388			NA	NA	NA	NA				NA	
										NA	NA	NA	NA				NA	
Healthcare Organizations	AGGREGATE DISCLOSURE																	
	Aggregate amount attributable to transfers of value to such Recipients				Only Third Parties		NA	NA	NA	NA	NA	NA	NA	NA	NA			
	Number of Recipients in aggregate disclosure				Only Third Parties		NA	NA	NA	NA	NA	NA	NA	NA	NA			
	% of the number of Recipients included in the aggregate disclosure in the total number of Recipients disclosed				Only Third Parties		NA	NA	NA	NA	NA	NA	NA	NA	NA			

* In the case of direct support for the formation of HCPs that they exercise in a private context, it will be necessary to publish the data in individual or aggregate form, depending on whether the HCP has given consent