

PUBBLICAZIONE 2021 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)																
	Soverini Simona	Bologna	Italy	Dipartimento di Medicina Specialistica, Diagnostica e	NA	NA	NA*	NA*	NA*					800		NA	800
	Marchini Sergio	Rozzano	Italy	Campus Building C. Via Manzoni 56, 20089 Rozzano	NA	NA	NA*	NA*	NA*					640		NA	640
					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*					2,450	1,089	NA	3,539
	Number of Recipients in aggregate disclosure				NA	NA	NA*	NA*	NA*							NA	2
	% of the number of Recipients included in the aggregate disclosure in the total number of Recipients disclosed				NA	NA	NA*	NA*	NA*							NA	50%
Healthcare Third Parties	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)																
	Selene S.r.l. – Eventi e Congressi	Torino	Italy	Via G. MEDICI 23, 10143 Torino	Only HCO	NA	10,440			NA	NA	NA	NA			NA	10,440
	SIMTIIPRO Srl	Milano	Italy	Via desiderio 21, 20131 Milano	Only HCO	NA	10,750			NA	NA	NA	NA			NA	10,750
	M2 Congressi srl	Milano	Italy	Via Carlo Farini, 81, 20159 Milano MI	Only HCO	NA	25,728			NA	NA	NA	NA			NA	25,728
	LC CONGRESSI S.R.L	Padova	Italy	VIA EUGANEA 45, 35141 PADOVA, PD, IT	Only HCO	NA	2,000			NA	NA	NA	NA			NA	2,000
	Scuola Medica Ospedaliera	Roma	Italy	Borgo Santo Spirito ,3 00193 Roma (RM)	Only HCO	NA	1,000			NA	NA	NA	NA			NA	1,000
	OIC GROUP	Firenze	Italy	Viale Giacomo Matteotti, 7, 50121 Firenze FI	Only HCO	NA	3,500			NA	NA	NA	NA			NA	3,500
										NA	NA	NA	NA			NA	
										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